

Diabetes

Insulin Signaling

PTPases

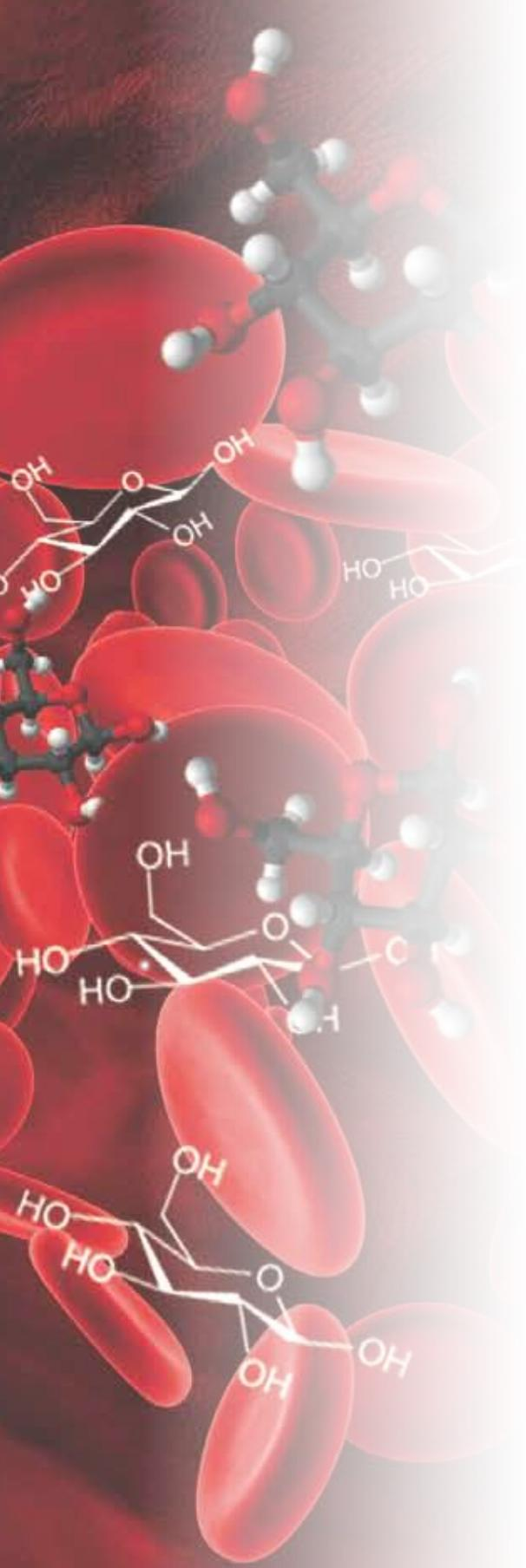
PI 3-K / Akt Pathway

GSK-3

Nutrient Sensing (mTOR / AMPK)

PPARs

GLP-1, DPPIV & Neprilysin



Enabling Discovery in Life Science®

ENZO LIFE SCIENCES, INC.

Enzo Life Sciences, Inc., a subsidiary of Enzo Biochem, Inc., is organized to lead in the development, production, marketing, and sale of innovative life science research reagents worldwide. Now incorporating the skills, experience, and products of ALEXIS Biochemicals, acquired in 2007, BIOMOL International, acquired in 2008, and ASSAY DESIGNS, acquired in 2009, Enzo Life Sciences provides over 25 years of business experience in the supply of research biochemicals, assay systems and biological reagents “Enabling Discovery in Life Science®”.

Based on a very substantial intellectual property portfolio, Enzo Life Sciences, Inc. is a major developer and provider of advanced assay technologies across research and diagnostic markets. A strong portfolio of labeling probes and dyes provides life science environments with tools for target identification and validation, and high content analysis via gene expression analysis, nucleic acid detection, protein biochemistry and detection, molecular biology, and cellular analysis.

- **Genomic Analysis**
- **Cellular Analysis**
- **Post-translational Modification**
- **Signal Transduction**
- **Cancer & Immunology**
- **Drug Discovery**

In addition to our wide range of catalog products, a complementary range of highly specialized custom services are also offered to provide tailor-made solutions for researchers. These include small molecule organic synthesis, peptide synthesis, protein expression, antibody production and immunoassay development.

International Edition 06/2010

www.enzolifesciences.com

Content

Introduction	4-5
Insulin Signaling	6-9
Protein Tyrosine Phosphatases [PTPases]	10-11
PI 3-Kinase [PI 3-K] Pathway	12-13
Akt [PKB; Protein Kinase B]	14-17
GSK-3 [Glycogen Synthase Kinase-3]	18-21
Nutrients & Nutrient Sensing mTOR Pathway AMPK Pathway	22-25
PPARs [Peroxisome Proliferator-activated Receptor]	26-29
Proteolytic Regulation of GLP-1 by DPPIV and Neprilysin	30-32
Antidiabetic Agents	33
HDACs	34-35
Sirtuins [SIRT6]	36-37
International Distributors	38-39

Introduction

Insulin and Diabetes

Diabetes is a chronic metabolic disorder affecting ~5% of the population in industrialized nations. The disease can be divided into two major types referred to as type 1 and type 2. Type 1 diabetes results from autoimmune destruction of pancreatic β -cells. This leads to a complete dependence on exogenous insulin to regulate blood glucose levels. The presence of auto-islet antibodies (glutamic acid decarboxylase (GAD), insulin and insulinoma antigen-3) is the best current marker to distinguish between diabetes type 1 and type 2.

The more prevalent form, type 2 diabetes, accounts for more than 90% of patients. Type 2 diabetes is characterized by the abnormal metabolism of glucose and fat. Several studies have established that type 2 diabetes is caused by a combination of insulin resistance in skeletal muscle, liver and adipose tissue and impaired insulin secretion from the pancreatic β -cells. Pancreatic β -cell dysfunction is a key element in the pathogenesis of type 2 diabetes. The reasons for β -cell failure are not completely understood. Several theories exist to explain the decrease in insulin secretion. These include a reduction of β -cell mass and impaired β -cell function through either β -cell exhaustion, desensitization to glucose stimulation, deposition of islet amyloid or lipotoxicity. With time, progressive β -cell failure leads to insulin deficiency and hyperglycemia. Hyperglycemia is the hallmark of type 2 diabetes and contributes to the pathogenesis of diabetes by impairing both insulin action and insulin secretion. Further progression of diabetes leads to the development of diabetic complications which include various microvascular (retinopathy, nephropathy and neuropathy) and macrovascular (cardiovascular and cerebrovascular) disorders.

A model for the development of type 2 diabetes was proposed by J.M. Fernandez-Real & J.C. Pickup (Figure 1). In this model several environmental risk factors such as diet, obesity, exercise, stress and age are thought to reduce insulin release and action in individuals with a susceptibility to the disease.

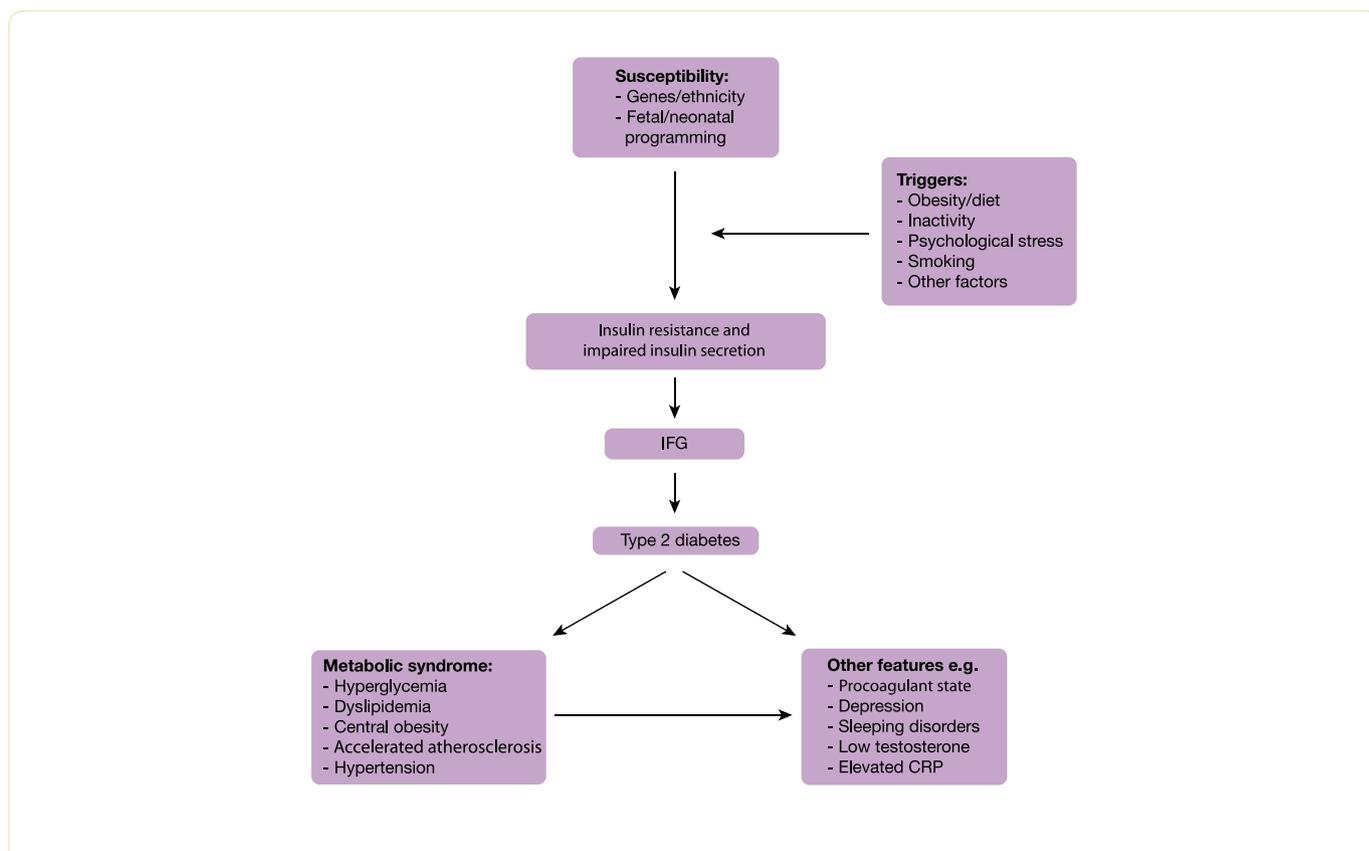


FIGURE 1: A model for the development of type 2 diabetes. Diabetes and the metabolic syndrome develop when environmental trigger factors, such as obesity, cause insulin resistance and impaired insulin secretion in genetically or otherwise susceptible people. Abbreviation: IGT/IFG, impaired glucose tolerance/impaired fasting glucose. Adapted from: Innate immunity, insulin resistance and type 2 diabetes: J.M. Fernandez-Real & J.C. Pickup; Trends Endocrinol. Metab. **19**, 10 (2008).

In addition it was suggested that a disorder of the innate immune system is responsible for the pathophysiology of type 2 diabetes, insulin resistance and several other chronic diseases such as atherosclerosis. In the proposed model (Figure 2), type 2 diabetes is associated with a general activation of the innate immune system, in which there is a chronic, cytokine-mediated state of low grade inflammation. This would explain most pathophysiologies of type 2 diabetes, including insulin resistance and impaired β -cell function, and many of the associated clinical conditions such as atherosclerosis and dyslipidemia.

LITERATURE REFERENCES:

The twentieth century struggle to decipher insulin Signaling: P. Cohen; Nat. Rev. Mol. Cell Biol. **7**, 867 (2006) (Review) ■ Role of insulin, adipocyte hormones, and nutrient-sensing pathways in regulating fuel metabolism and energy homeostasis: a nutritional perspective of diabetes, obesity, and cancer: S. Marshall; Sci. STKE **2006**, re7 (2006) (Review) ■ Abnormalities in insulin secretion in type 2 diabetes mellitus: P.J. Guillausseau, et al.; Diabetes Metab. **34 Suppl 2**, S43 (2008) (Review) ■ Beta-cell failure in type 2 diabetes mellitus: C. Lencioni, et al.; Curr. Diab. Rep. **8**, 179 (2008) (Review) ■ Genes and type 2 diabetes mellitus: L. Groop & V. Lyssenko; Curr. Diab. Rep. **8**, 192 (2008) (Review) ■ Innate immunity, insulin resistance and type 2 diabetes: J.M. Fernandez-Real & J.C. Pickup; Trends Endocrinol. Metab. **19**, 10 (2008) ■ Islet inflammation in type 2 diabetes: from metabolic stress to therapy: M.Y. Donath, et al.; Diabetes Care **31 Suppl. 2**, S161 (2008) ■ Inflammatory mechanisms in the regulation of insulin resistance: H. Tilg & A.R. Moschen; Mol. Med. **14**, 222 (2008)

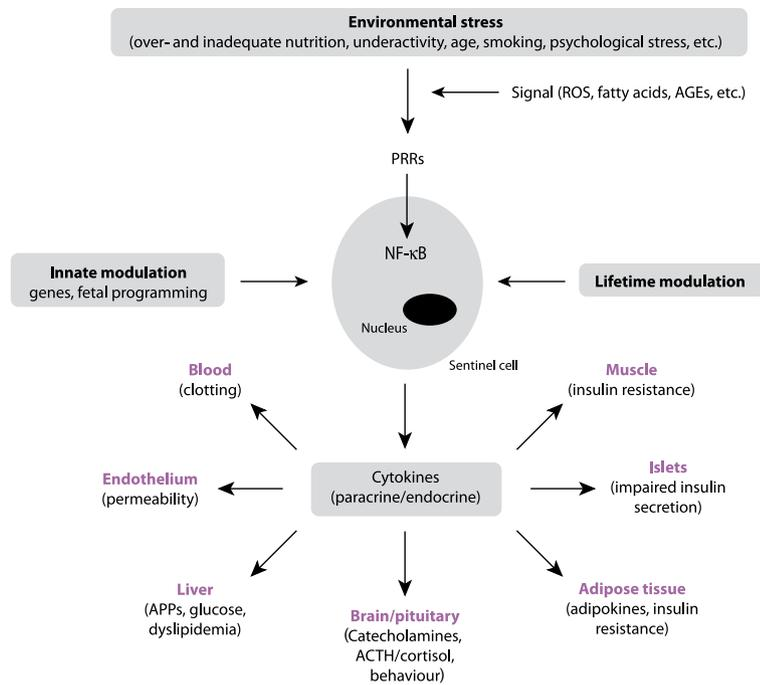


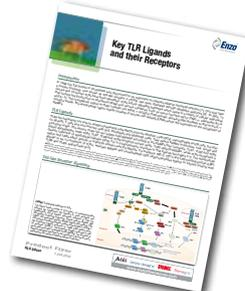
FIGURE 2: Innate immunity and type 2 diabetes. Sentinel cells, such as macrophages, hepatocytes, endothelium and adipocytes, detect environmental threats to the host through pattern-recognition receptors (PRRs). This activates the inflammatory NF- κ B pathway, releasing cytokines that act on many cells in the body to produce the clinical and biochemical features of the metabolic syndrome and type 2 diabetes (in parentheses). Abbreviations: ACTH, adrenocorticotrophic hormone; AGE, advanced glycosylation end product; APP, acute-phase protein; ROS, reactive oxygen species. Adapted from: Innate immunity, insulin resistance and type 2 diabetes: J.M. Fernandez-Real & J.C. Pickup; Trends Endocrinol. Metab. **19**, 10 (2008).

Latest Insight

Recent findings indicate that TLRs (Toll-like receptors) - especially TLR4, a sensor for endogenous lipids and free fatty acids - mediate the link between the metabolic and immune systems. Studies on TLR4 deficient mice have shown that these animals become obese, apparently due to overeating, demonstrating once more the important role of TLR4.

LITERATURE REFERENCES:

Toll-like receptors in endocrine disease and diabetes: W. Kanczkowski, et al.; Neuroimmunomodulation **15**, 54 (2008)



Visit www.enzolifesciences.com for a complete listing or ask for a free copy of our product flyers.

Insulin Signaling

The anabolic hormone insulin is released by the β -cells of the pancreas islets in response to elevated levels of nutrients in the blood. It serves as the key regulator of plasma glucose by maintaining a balance between the intestinal absorption, liver production and cellular uptake of glucose and its metabolism by peripheral tissues. Insulin triggers the uptake of glucose, fatty acids and amino acids into adipose tissue, muscle and the liver and promotes the storage of these nutrients in the form of glycogen, lipids and proteins. At times of starvation, insulin levels are low and this directs insulin-sensitive cells to mobilize stored fuels.

At a cellular level, insulin/insulin-like growth factor (IGF)-signaling regulates multiple processes, including carbohydrate and lipid metabolism, gene transcription, DNA synthesis, apoptosis, and cell proliferation. Insulin and insulin-like growth factors (IGF-1 and IGF-2) bind to members of the insulin receptor tyrosine kinase family, including insulin receptors (IRs) and the IGF-1 receptor (IGF1R). The heterotetrameric insulin receptor consists of 2 extracellular β subunits that bind insulin and 2 transmembrane α subunits with tyrosine kinase activity. Upon ligand binding, the activated insulin receptor phosphorylates tyrosine residues on intracellular substrates such as the insulin receptor substrates (IRS1-4), IRS5/DOK4, IRS/DOK5, Gab-1 and Shc isoforms. This leads to the activation of downstream signaling molecules including the mitogen-activated protein kinase (MAPK) cascade, other serine/threonine kinases (Akt, GSK-3, mTOR, PKC) and phosphoinositide 3-kinase (PI 3-K).

LITERATURE REFERENCE:

Insulin signaling and the regulation of glucose transport: L. Chang, et al.; Mol. Med. **10**, 65 (2004) • Insulin signaling: P. Bevan; J. Cell Sci. **114**, 1429 (2001)

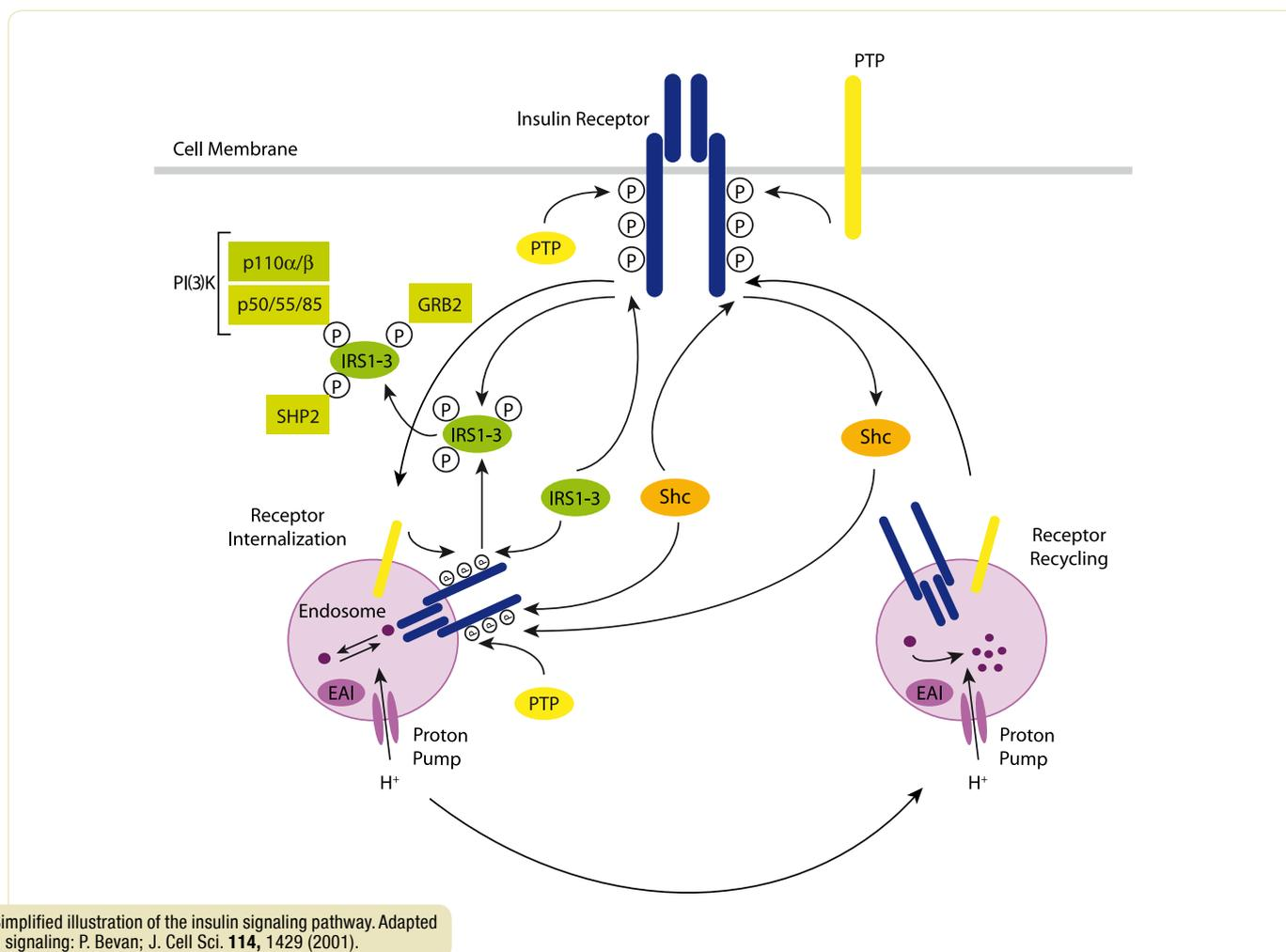


FIGURE 3: Simplified illustration of the insulin signaling pathway. Adapted from: Insulin signaling: P. Bevan; J. Cell Sci. **114**, 1429 (2001).

Insulin

Insulin, mAb (SPM139)

ADI-905-433-1

1 ml

CLONE: SPM139. ISOTYPE: Mouse IgG1. IMMUNOGEN: Swine insulin. SPECIFICITY: Recognizes human, rat, bovine, rabbit and pig Insulin. APPLICATION: IHC.

C-Peptide (human), mAb (91/1F9)

ALX-803-049-C100

100 µg

CLONE: 91/1F9. ISOTYPE: Mouse IgG2b. IMMUNOGEN: Human proinsulin. SPECIFICITY: Recognizes human C-peptide. Cross-reacts with human proinsulin. APPLICATION: IHC (FS), IP, WB.

Insulin Receptor (IR) Peptides

Insulin receptor (1142-1153)

[H-Thr-Arg-Asp-Ile-Tyr-Glu-Thr-Asp-Tyr-Tyr-Arg-Lys-OH]

BML-P314-0500

0.5 mg

This sequence, is derived from the insulin receptor β -subunit cytoplasmic domain, inclusive of regulatory autophosphorylation sites Tyr^{1146/1150/1151}. It has been used as an insulin receptor tyrosine kinase substrate and the phosphorylated peptide as a protein tyrosine phosphatase substrate.

[pTyr¹¹⁴⁶]Insulin receptor (1142-1153)

[H-Thr-Arg-Asp-Ile-pTyr-Glu-Thr-Asp-Tyr-Tyr-Arg-Lys-OH]

BML-P315-0500

0.5 mg

This phosphopeptide sequence, corresponds to the insulin receptor regulatory region. Can be used as a tyrosine phosphatase substrate.

[pTyr^{1146/1150}]Insulin receptor (1142-1153)

[H-Thr-Arg-Asp-Ile-pTyr-Glu-Thr-Asp-pTyr-Tyr-Arg-Lys-OH]

BML-P321-0500

0.5 mg

This sequence, is derived from the insulin receptor β -subunit cytoplasmic domain, inclusive of regulatory autophosphorylation sites Tyr^{1146/1150/1151}. The unphosphorylated version of this peptide (Prod. Nr. BML-P314) has been used as an insulin receptor kinase substrate. The phosphorylated peptide has been used as a protein tyrosine phosphatase substrate.

[pTyr^{1146/1150/1151}]Insulin receptor (1142-1153)

[H-Thr-Arg-Asp-Ile-pTyr-Glu-Thr-Asp-pTyr-pTyr-Arg-Lys-OH]

BML-P318-0500

0.5 mg

This phosphopeptide can be used as a substrate for protein tyrosine phosphatases.

[pTyr¹¹⁵⁰]Insulin receptor (1142-1153)

[H-Thr-Arg-Asp-Ile-Tyr-Glu-Thr-Asp-pTyr-Tyr-Arg-Lys-OH]

BML-P316-0500

0.5 mg

This phosphopeptide corresponds to the insulin receptor regulatory region. Can be used as a tyrosine phosphatase substrate.

[pTyr¹¹⁵¹]Insulin receptor (1142-1153)

[H-Thr-Arg-Asp-Ile-Tyr-Glu-Thr-Asp-Tyr-pTyr-Arg-Lys-OH]

BML-P317-0500

0.5 mg

This phosphopeptide corresponds to the insulin receptor regulatory region. Can be used as a tyrosine phosphatase substrate.

Insulin receptor phosphopeptide, sampler pack

BML-P319-0001

1 Kit

Contains 0.5 mg each of the insulin receptor peptides (BML-P314, BML-P315, BML-P316, BML-P317, BML-P318 and BML-P321).

Insulin Receptor (IR) Antibodies

Product	Specificity	Application	Prod. No.	Size
[pTyr ¹¹⁵⁰ /Tyr ¹¹⁵¹]Insulin receptor, mAb (10C3)	Human	ELISA, WB	ADI-905-645-100	100 µg
[pTyr ^{1158/1162/1163}]Insulin receptor and [pTyr ^{1131/1135/1136}]IGF-1 receptor, pAb	Human	WB	BML-SA392-0020 BML-SA392-0100	20 µl 100 µl
[pTyr ¹¹⁵⁸]Insulin receptor and [pTyr ¹¹³¹]IGF-1 Receptor (human), pAb	Human	WB	BML-SA390-0020 BML-SA390-0100	20 µl 100 µl
[pTyr ¹³²²]Insulin receptor, mAb (21G12)	Human, mouse and dog	ELISA, WB	ADI-905-646-100	100 µg
[pTyr ⁹⁷²]Insulin receptor, pAb	Human, mouse and (rat)	WB	ADI-CSA-720-E	100 µl
Insulin receptor β subunit, mAb (CT-3)	Human, mouse, rat and monkey	IHC, WB	BML-SA432-0100	100 µg
Insulin receptor β subunit, mAb (C18C4)	Human	WB	ADI-905-683-100	100 µg
Insulin receptor, mAb (9H4)	Human, mouse, rat and dog	ELISA, IHC, WB	ADI-905-647-100	100 µg

Insulin Receptor Substrates (IRS) Peptide and Antibodies

IRS-1 (Tyr⁶⁰⁸) peptide

[H-Lys-Lys-His-Thr-Asp-Asp-Gly-Tyr-Met-Pro-Met-Ser-Pro-Gly-Val-Ala-OH]

BML-P320-0001 1 mg

This sequence, is derived from (insulin receptor substrate-1) (IRS-1) inclusive of Tyr-608 (mouse) -612 (human). It contains the insulin receptor tyrosine kinase substrate motif YMXM (Tyr-Met-X-Met). This peptide has been used as a substrate for purified insulin receptor ($K_m=90 \mu\text{M}$) and other tyrosine kinases in phosphocellulose binding assays. The tyrosine phosphorylated version of this peptide binds to phosphatidylinositol 3-kinase (PI 3-K) SH2 domain and activates the enzyme.

[pTyr⁶¹²]IRS-1, pAb

ADI-CSA-725-E 100 μl

From rabbit. **IMMUNOGEN:** Synthetic phospho-peptide derived from the sequence of human IRS-1 (sequence identical with mouse, rat and chicken). **SPECIFICITY:** Recognizes human, (mouse, rat and chicken) IRS-1. **APPLICATION:** WB.

[pTyr¹²²⁹]IRS-1 (human), pAb

BML-SA393-0020 20 μl

BML-SA393-0100 100 μl

From rabbit. **IMMUNOGEN:** Synthetic peptide corresponding to a portion of human IRS-1 (insulin receptor substrate-1) phosphorylated at Tyr¹²²⁹. **SPECIFICITY:** Recognizes human IRS-1 phosphorylated at Tyr¹²²⁹. **APPLICATION:** WB. **BP:** BML-SP393.

[pTyr⁹⁴¹]IRS-1 (human), pAb

BML-SA394-0020 20 μl

BML-SA394-0100 100 μl

From rabbit. **IMMUNOGEN:** Synthetic peptide corresponding to a portion of human IRS-1 (insulin receptor substrate-1) phosphorylated at Tyr⁹⁴¹. **SPECIFICITY:** Recognizes human IRS-1 phosphorylated at Tyr⁹⁴¹. Tyrosine 941 of IRS-1 is in a phosphatidylinositol 3-kinase (PI 3-K) binding site motif. **APPLICATION:** WB. **BP:** BML-SP394.

[pTyr⁸⁹⁶]IRS-1 (human), pAb

BML-SA396-0020 20 μl

BML-SA396-0100 100 μl

From rabbit. **IMMUNOGEN:** Synthetic peptide corresponding to a portion of human IRS-1 (insulin receptor substrate-1) phosphorylated at Tyr⁸⁹⁶. **SPECIFICITY:** Recognizes human IRS-1 phosphorylated at Tyr⁸⁹⁶. Tyrosine 896 of IRS-1 is a Grb-2 SH2 domain binding site. **APPLICATION:** WB. **BP:** BML-SP396.

Insulin Receptor (IR) Inhibitors

Hydroxy-2-naphthalenylmethylphosphonic acid

[HNMPA]

BML-EI247-0005 5 mg

BML-EI247-0025 25 mg

Membrane impermeable inhibitor of insulin receptor tyrosine kinase activity ($IC_{50}=100\mu\text{M}$). Was shown to inhibit both tyrosine and serine autophosphorylation by the human insulin receptor.

HNMPA-(AM)₃

BML-EI248-0005 5 mg

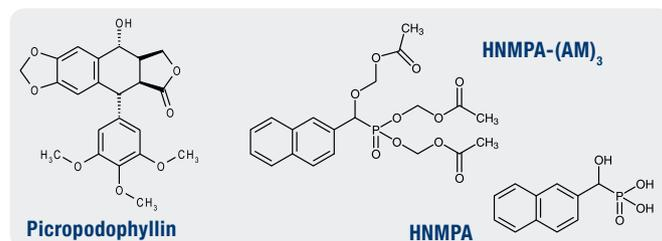
BML-EI248-0025 25 mg

Cell permeable analog of HNMPA. Inhibits insulin receptor tyrosine kinase activity ($IC_{50}=100\mu\text{M}$) and insulin-stimulated glucose oxidation in isolated rat adipocytes ($IC_{50}=10\mu\text{M}$). Has no effect on PKA (at concs. up to 1mM) or PKC (at concs. up to 420 μM).

Picropodophyllin

BML-EI372-0001 1 mg

Potent and selective inhibitor of insulin-like growth factor 1 receptor (IGF-1R, $IC_{50}=6\text{nM}$). Efficiently blocks IGF-1R activity, reduces pAKT and pERK1/2, induces apoptosis in IGF-1R-positive tumor cells and causes complete tumor regression in xenografted and allografted mice. Picropodophyllin downregulates IGF-1R by interfering with the action of β -arrestin 1/MDM2.



Insulin Receptor (IR) Kinase

Insulin receptor kinase β subunit (human), (rec.) (GST-tag)

BML-SE195-0020 20 μg

Produced in insect cells. The 46kDa cytoplasmic domain of the β -subunit of the human insulin receptor (βIRK) (aa 941-1343) is fused to a GST-tag. **APPLICATION:** Enzyme kinetics and inhibition studies and for phosphorylation of target substrates.

Insulin receptor kinase β subunit (human) (rec.) (His-tag)

BML-SE516-0020 20 μg

Produced in insect cells. The cytoplasmic (kinase) domain of human insulin receptor is fused to a His-tag.

Shc

Shc, pAb

ALX-210-535-R100 100 µl

From rabbit. **IMMUNOGEN:** Recombinant human Shc portion (aa 1180-1620). **SPECIFICITY:** Recognizes human, mouse and hamster p46shc, p52shc and p66shc isoforms of phosphorylated and unphosphorylated Shc. **APPLICATION:** ELISA, IP, WB.

Shc, mAb (11F6)

ADI-905-665-100 100 µg

CLONE: 11F6. **ISOTYPE:** Mouse IgG1. **IMMUNOGEN:** Synthetic peptide derived from sequence near the carboxy-terminus of Shc, conjugated to hemocyanin. **SPECIFICITY:** Recognizes human and mouse Shc. **APPLICATION:** ICC, IHC, WB.

[pSer³⁶]Shc, mAb (6E10)

ALX-804-358-C100 100 µg

CLONE: 6E10. **ISOTYPE:** Mouse IgG1. **IMMUNOGEN:** Synthetic peptide corresponding to aa 33-39 (E³³LPpSPSA³⁹) of shc phosphorylated at Ser³⁶. **SPECIFICITY:** Recognizes human, mouse and dog shc phosphorylated at Ser³⁶. Does not cross-react with non-phosphorylated shc. **APPLICATION:** WB.

[pTyr³¹⁷]Shc, mAb (15E11)

ALX-804-359-C100 100 µg

CLONE: 15E11. **ISOTYPE:** Mouse IgG1. **IMMUNOGEN:** Synthetic peptide corresponding to a portion of Shc phosphorylated at Tyr³¹⁷. **SPECIFICITY:** Recognizes human, mouse and dog Shc phosphorylated at Tyr³¹⁷. **APPLICATION:** IHC, WB.

[pTyr^{239/240}]Shc, mAb (1E3)

ALX-804-357-C100 100 µg

CLONE: 1E3. **ISOTYPE:** Mouse IgG1. **IMMUNOGEN:** Synthetic peptide corresponding to aa 236-243 (D²³⁶HQpYpYND²⁴³) of shc phosphorylated at Tyr²³⁹ and Tyr²⁴⁰. **SPECIFICITY:** Recognizes human, mouse and dog p46shc, p52shc and p66shc when phosphorylated at Tyr^{239/240}. Does not cross-react with non-phosphorylated shc. **APPLICATION:** IHC, WB.

Insulin-like Growth Factor

Insulin-like growth factor (IGF)-1 is a powerful mitogenic peptide synthesized predominantly by the liver. It circulates in the plasma and is one of the main mediators of the actions of growth hormone (GH). Postnatal, IGF-1 promotes cell proliferation, differentiation and survival. However, IGF-1 has also effects on insulin sensitivity and several studies suggest that IGF-1 levels may be altered in insulin resistance states and type 2 diabetes.

LITERATURE REFERENCE:

The role of IGF-1 and its binding proteins in the development of type 2 diabetes and cardiovascular disease: V.A. Ezzat, et al.; *Diabetes Obes. Metab.* **10**, 198 (2008) (Review)

IGF-1 (human), (rec.)

ADI-908-059-020 20 µg

Recombinant human IGF-1 protein expressed in *E. coli*.

[pTyr¹³¹⁶]IGF-1 receptor, mAb (2B9)

ADI-905-643-100 100 µg

CLONE: 2B9. **ISOTYPE:** Mouse IgG1. **IMMUNOGEN:** Synthetic phospho-peptide derived from the sequence of human IGF-1 receptor, conjugated to hemocyanin. **SPECIFICITY:** Recognizes human IGF-1 receptor. **APPLICATION:** ELISA, WB.

IGF-1 receptor kinase β (human), (rec.) (GST-tag)

BML-SE232-0020 20 µg

Produced in baculovirus. The 46kDa cytoplasmic domain of the β-subunit of human IGF-1 receptor kinase (IGF-1RK) (aa 929-1337) is fused to GST. **APPLICATION:** Kinetic studies, inhibitor screening, and phosphorylation of target substrates.

IGF-1 receptor, mAb (7G11)

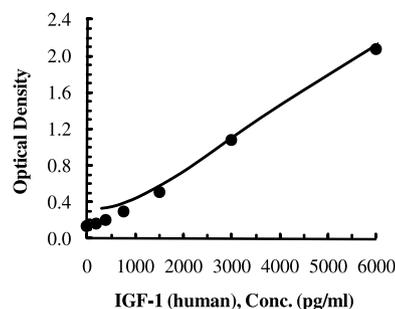
ADI-905-644-100 100 µg

CLONE: 7G11. **ISOTYPE:** Mouse IgG1. **IMMUNOGEN:** Synthetic peptide derived from the sequence near the carboxy-terminus of human IGF-1 receptor, conjugated to hemocyanin. **SPECIFICITY:** Recognizes human IGF-1 receptor. **APPLICATION:** WB.

IGF-1 (human), EIA kit

ADI-900-150 96 wells

For the quantitative determination of human IGF-1 in plasma and serum. **SENSITIVITY:** 48.5 pg/ml (range 187 - 6,000 pg/ml).



HIGHLIGHT

Protein Tyrosine Phosphatases [PTPases]

Protein tyrosine phosphatases (PTPases) are important in cellular signal transduction and metabolism. The PTPase family can be divided into two categories: Intracellular (non-receptor type) PTPases (e.g. PTP1B); and transmembrane (receptor type) PTPases (e.g. LAR). Several PTPases have been described to catalyze the dephosphorylation of the insulin receptor. These include PTP α , LAR, CD45, PTP ϵ , SHP2, TC-PTP and PTP1B. However, the implication of these PTPases in the regulation of insulin signaling is still controversial. The most convincing data supports a critical role for PTP1B in insulin action, which makes PTP1B an attractive candidate for the drug design for type 2 diabetes and obesity.

LITERATURE REFERENCES:

Protein tyrosine phosphatases: the quest for negative regulators of insulin action: E. Asante-Appiah & B.P. Kennedy; Am. J. Physiol. Endocrinol. Metab. **284**, E663 (2003) (Review) • Involvement of the small protein tyrosine phosphatases TC-PTP and PTP1B in signal transduction and diseases: from diabetes, obesity to cell cycle, and cancer: N. Dube & M.L. Tremblay; Biochim. Biophys. Acta **1754**, 108 (2005) • Protein tyrosine phosphatases: structure-function relationships: L. Taberner, et al.; FEBS J. **275**, 867 (2008) (Review)

PTP1B

PTP1B (human), (rec.)

BML-SE332-0050 50 μ g

Produced in *E. coli*. Human PTP1B (protein tyrosine phosphatase 1B) (aa 1-322). **APPLICATION:** Useful for the study of tyrosine phosphatase kinetics, substrate specificity and for screening inhibitors.

[pTyr⁹⁹²]EGF receptor (988-998)

[EGFR (988-998); H-Asp-Ala-Asp-Glu-pTyr-Leu-Ile-Pro-Gln-Gln-Gly-OH]

BML-P323-0001 1 mg

This sequence, DADEpYLIPQQG, is from an autophosphorylation site (Tyr⁹⁹²) of epidermal growth factor receptor (EGFR). An excellent substrate for mammalian PTP1B with $K_m=3.9 \mu$ M) and *Yersinia* PTPase. May be used in non-radioactive phosphatase assays incorporating the BIOMOL GREEN™ reagent (Prod. Nr. BML-AK111).

PTP1B drug discovery kit

BML-AK822-0001 1 Kit

The PTP1B Tyrosine Phosphatase Drug Discovery Kit is a colorimetric, non-radioactive assay designed to measure the phosphatase activity of purified PTP1B. This 96 well assay is useful for screening inhibitors and modulators of PTP1B activity. The kit includes human, recombinant PTP1B (residues 1-322; MW=37.4 kDa), expressed in *E. coli*. The detection of free-phosphate released is based on the classic Malachite green assay and offers the advantages of convenient, 1-step detection and excellent sensitivity, without radioactivity.

LAR

LAR (human), (rec.)

BML-SE113-0200 200 U

Produced in *E. coli*. Soluble catalytic LAR-D1 domain (350 aa).

LAR Tyrosine Phosphatase Assay Kit for Drug Discovery

BML-AK815-0001 1 Kit

The LAR tyrosine phosphatase assay kit is a colorimetric, non-radioactive assay designed in a 96 well microplate format. It may be used to screen inhibitors and modulators of LAR phosphatase activity, and perform other experiments on LAR enzyme kinetics. The kit includes purified human LAR cytoplasmic domain (residues 1275-1613; MW=37 kDa), expressed in an *E. coli* expression system. The assay entails LAR dephosphorylation of a synthetic 12-mer peptide based on the regulatory autophosphorylation region of the IR β subunit cytoplasmic domain (residues 1142-1153) and phosphorylated at LAR's preferred dephosphorylation site Tyr-1146. The phosphate released is quantified with the BIOMOL GREEN™ reagent. A tyrosine phosphatase inhibitor, RWJ-60475 is included as a control "hit" for inhibitor screening.



HIGHLIGHT

TC-PTP

T cell protein tyrosine phosphatase (human), (rec.)

BML-SE114-0200 200 U
Produced in *E. coli*. T cell phosphatase with a C-terminal 11 kDa deletion (TCΔC11) (aa 1-317).

PTP-IA2 [Protein Tyrosine Phosphatase IA2]

PTP-IA2 (catalytic domain) (human), (rec.) (GST-tag)

BML-SE278-0020 20 µg
Produced in *E. coli*. The catalytic domain of human PTP-IA2 (protein tyrosine phosphatase IA2) (aa 693-970) is fused at the N-terminus to a GST-tag.
APPLICATION: Study of enzyme kinetics and regulation, dephosphorylation of target substrates, and inhibitor screening.

PTP-IA2 (intracellular domain) (618-628), mAb (98/4H6)

ALX-803-056-C100 100 µg
CLONE: 98/4H6. **ISOTYPE:** Mouse IgG1. **IMMUNOGEN:** Recombinant human IA-2ic (protein tyrosine phosphatase IA2 (intracellular domain)). **SPECIFICITY:** Recognizes an epitope (aa 618-628) of human, mouse and rat IA-2ic. **APPLICATION:** IHC (FS), IP, WB.

PTP-IA2 (intracellular domain) (605-772), mAb (97/4D9)

ALX-803-057-C100 100 µg
CLONE: 97/4D9. **ISOTYPE:** Mouse IgG1. **IMMUNOGEN:** Recombinant human IA-2ic (protein tyrosine phosphatase IA2 (intracellular domain)). **SPECIFICITY:** Recognizes an epitope (aa 605-772) of human, mouse and rat IA2ic. **APPLICATION:** IHC (FS), IP, WB.

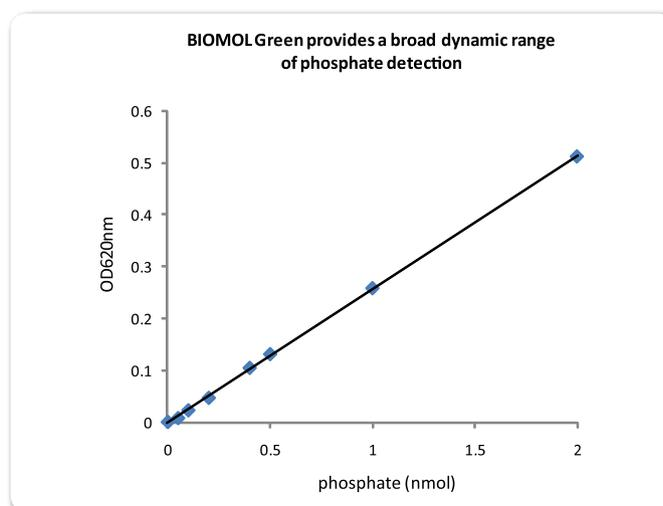
PTP-IA2 (intracellular domain) (771-979), mAb (103/2C4)

ALX-803-058-C100 100 µg
CLONE: 103/2C4. **ISOTYPE:** Mouse IgG1. **IMMUNOGEN:** Recombinant human IA-2ic (protein tyrosine phosphatase IA2 (intracellular domain)). **SPECIFICITY:** Recognizes an epitope (aa 771-979) of human, mouse and rat IA2ic. **APPLICATION:** IHC (FS), IP, WB.

Related Products

BIOMOL Green™ Reagent

BML-AK111-0250 250 ml
BIOMOL GREEN™ provides a simple and convenient method for colorimetric phosphate quantitation (abs. 600-680 nm). Unlike other molybdate/malachite green-based assays, BIOMOL GREEN™ doesn't require multiple solutions or reagents prepared fresh on the day of the assay. BIOMOL GREEN™ is extremely stable (>6 mos. at 4°C) and is simply mixed, at room temperature, with any enzymatic reaction that has released free phosphate. Reported applications include assays for phospholipid phosphatases, tyrosyl-tRNA synthetase (coupled with pyrophosphatase) and a viral RNA triphosphatase. BIOMOL GREEN™ has perhaps been most widely applied to protein phosphatase assays. Typically, a protein phosphatase (e.g. PTP1B (Prod. No. BML-SE332) or LAR (Prod. No. BML-SE113)) is simply incubated with an appropriate phosphopeptide (e.g. IR5,9,10 (Prod. No. BML-P318)). Addition of BIOMOL GREEN™ stops the reaction and begins color development, which is read 20-30 min later. BIOMOL GREEN™ may be used in cuvette or microplate-based assays and is ideal for high-throughput applications. Each kit comes with a standardized phosphate solution for assay calibration.



PI 3-Kinase [PI 3-K] Pathway

The PI 3-K (phosphatidylinositol 3-kinase) pathway plays a pivotal role in the metabolic and mitogenic action of insulin. PI 3-K is a heterodimeric enzyme consisting of a p85 regulatory subunit as well as a p110 catalytic subunit. It gets activated by several hormones including growth factors and insulin. Upon activation, PI 3-K catalyzes the phosphorylation of phosphoinositols (e.g. the formation of PIP3 from PIP2), which subsequently leads to the recruitment of Akt (PKB; protein kinase B) to the cell membrane. Activated Akt regulates the activity of many downstream proteins involved in multiple aspects of cellular physiology (see Figure 4). Among others, Akt phosphorylates and regulates components of the glucose transporter 4 (GLUT4) complex, protein kinase C (PKC) isoforms and GSK-3, all of which are critical for insulin-mediated metabolic effects.

LITERATURE REFERENCES:

Phosphoinositide 3-kinase in disease: timing, location, and scaffolding: M.P. Wymann & R. Marone; *Curr. Opin. Cell Biol.* **17**, 141 (2005) (Review) • Targeting phosphoinositide 3-kinase: moving towards therapy: R. Marone, et al.; *Biochim. Biophys. Acta* **1784**, 159 (2008) (Review)

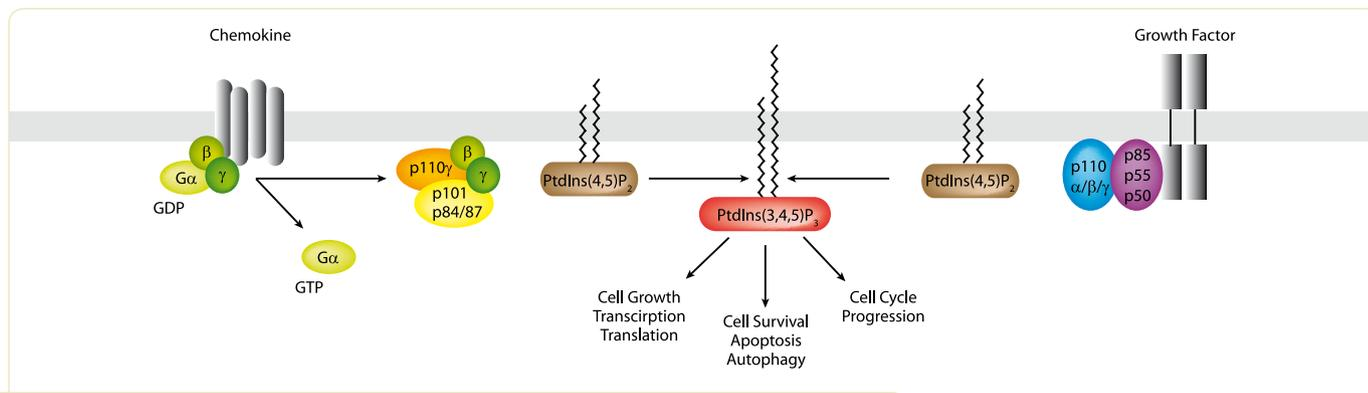


FIGURE 4: Simplified activation scheme of class I PI 3-Ks. Heterodimeric PI 3-K α , PI 3-K β , and PI 3-K δ complexes are activated downstream of growth factors, cytokine receptors and their substrates, whereas PI 3-K γ activation is triggered downstream of G protein-coupled receptors (GPCRs). Adapted from: *Targeting phosphoinositide 3-kinase: moving towards therapy*; R. Marone, et al.; *Biochim. Biophys. Acta* **1784**, 159 (2008).

PI 3-Kinase Proteins

PI 3-kinase p110 α /p85 α (human), (rec.)

BML-SE436-0020 20 μ g
Produced in insect cells. Complex of human full length p110 α and p85 α subunits of PI 3-kinase. **APPLICATION:** Kinetic and functional studies, phosphorylation of target substrates, drug screening.

PI 3-kinase p110 α /p85 (bovine), (rec.)

ALX-201-119-C005 5 μ g
Produced in Sf9 cells. Untagged bovine PI 3-K p110 α /p85 dimer.

PI 3-kinase p110 γ (human), (rec.)

ALX-201-055-C010 10 μ g
Produced in Sf9 cells. Full length human PI 3-kinase p110 γ fused at the N-terminus to a His-tag.

LIT: Cloning and characterization of a G protein-activated human phosphoinositide-3 kinase: B. Stoyanov, et al.; *Science* **269**, 690 (1995) • Lipid kinase and protein kinase activities of G-protein-coupled phosphoinositide 3-kinase gamma: structure-activity analysis and interactions with wortmannin: S. Stoyanova, et al.; *Biochem. J.* **324**, 489 (1997) • Linkage of G protein-coupled receptors to the MAPK signaling pathway through PI 3-kinase gamma: M. Lopez-Illasaca, et al.; *Science* **275**, 394 (1997) • Gbetagamma stimulates phosphoinositide 3-kinase-gamma by direct interaction with two domains of the catalytic p110 subunit: D. Leopoldt, et al.; *J. Biol. Chem.* **273**, 7024 (1998)

PI 3-Kinase Antibodies

PI 3-kinase, mAb (AB6)

ADI-KAM-PI200-E 100 μ g
CLONE: AB6. **ISOTYPE:** Mouse IgG1. **IMMUNOGEN:** Recombinant human p85 α protein. **SPECIFICITY:** Recognizes human, mouse and rat PI 3-kinase. **APPLICATION:** ICC, IP, WB.

PI 3-kinase p110 δ (human), pAb

ALX-210-747-C200 200 μ g
From rabbit. **IMMUNOGEN:** Synthetic peptide corresponding to aa 1026-1044 (S¹⁰²⁶WTKVNWLAHNVSKDNRQ¹⁰⁴⁴) of the C-terminal domain of human PI 3-kinase p110 δ . **SPECIFICITY:** Recognizes human PI 3-kinase p110 δ . Detects a band of ~110kDa by Western blot. **APPLICATION:** IHC (PS), ICC, WB. **BP:** ALX-165-034.

PI 3-Kinase Activators

PI 3-kinase SH3 domain binding peptide

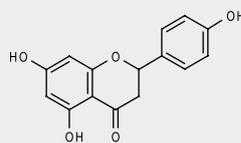
[H-Arg-Lys-Leu-Pro-Pro-Arg-Pro-Arg-Arg-OH]
BML-P311-0001 1 mg

IRS-1 (Tyr⁶⁰⁸) peptide

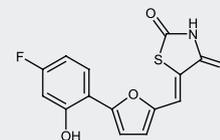
[H-Lys-Lys-His-Thr-Asp-Asp-Gly-Tyr-Met-Pro-Met-Ser-Pro-Gly-Val-Ala-OH]
BML-P320-0001 1 mg

PI 3-kinase Inhibitors

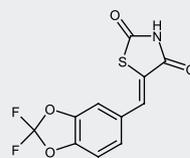
N-Acetyl-Asp-Tyr(2-malonyl)-Val-Pro-Met-Leu-NH₂ ALX-151-026-M001	1 mg
N-Acetyl-Asp-Tyr(PO₃H₂)-Val-Pro-Met-Leu-NH₂ ALX-151-027-M001	1 mg
(±)-Naringenin ALX-385-010-G001	1 g
AS-252424 ALX-270-465-M001	1 mg
ALX-270-465-M005	5 mg
ALX-270-465-M025	25 mg
AS-604850 ALX-270-461-M001	1 mg
ALX-270-461-M005	5 mg
ALX-270-461-M025	25 mg
AS-605240 ALX-270-462-M001	1 mg
ALX-270-462-M005	5 mg
ALX-270-462-M025	25 mg
Compound 15e ALX-270-455-M001	1 mg
ALX-270-455-M005	5 mg
Luteolin ALX-385-007-M010	10 mg
ALX-385-007-M050	50 mg
LY 294002 BML-ST420-0005	5 mg
BML-ST420-0025	25 mg
LY-303,511 ALX-270-410-M001	1 mg
ALX-270-410-M005	5 mg
PI-103 ALX-270-460-M001	1 mg
ALX-270-460-M005	5 mg
ALX-270-460-M025	25 mg
PI 3K-SH2-OMT BML-ST425-0500	0.5 mg
TGX-221 ALX-270-464-C100	100 µg
ALX-270-464-M001	1 mg
Wortmannin BML-ST415-0001	1 mg
BML-ST415-0005	5 mg
BML-ST415-0020	20 mg



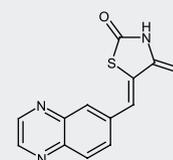
(±)-Naringenin



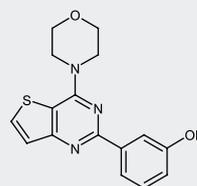
AS-252424



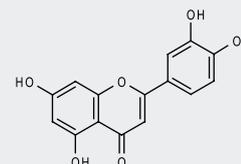
AS-604850



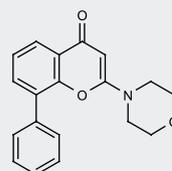
AS-605240



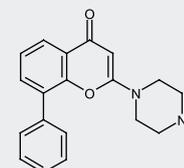
Compound 15e



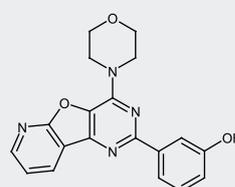
Luteolin



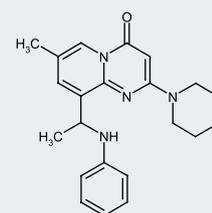
LY 294002



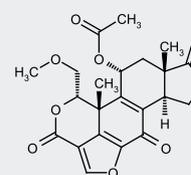
LY-303,511



PI-103



TGX-221



Wortmannin

Akt [PKB; Protein Kinase B]

Phosphorylation Site of the Akt Isoforms

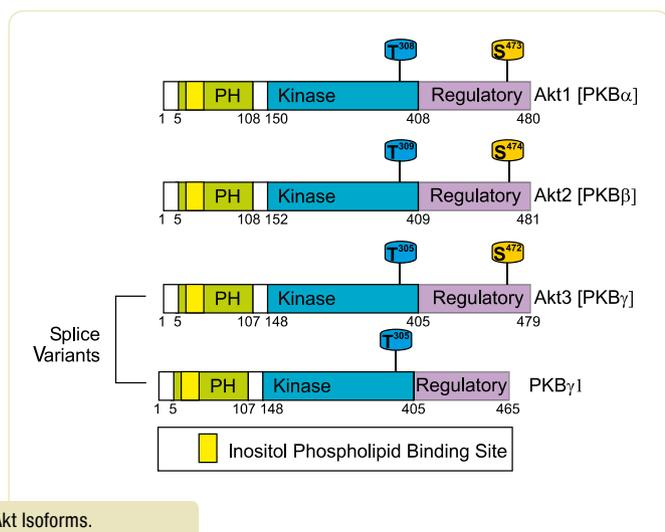


FIGURE 5: Akt Isoforms.

Akt1 shares 81% and 83% amino acid identity with Akt2 and Akt3, respectively. Although the three isoforms show broad tissue distribution, Akt1 is the most ubiquitously expressed. Akt2 is expressed at lower levels than Akt1 except in insulin-responsive tissues where it predominates. In particular, Akt2 is more abundant and more highly activated than Akt1 in adipocytes. Akt3 is expressed in testes and brain. The C-terminal sequence of the three isoforms are shown below. Note that the phosphorylation site differs slightly between the three isoforms. Thus the phosphorylation site Ser⁴⁷³ of Akt1 corresponds to Ser⁴⁷⁴ in Akt2 and Ser⁴⁷² in Akt3 respectively.

Akt1: V⁴⁶¹DSERRPHFPQFS⁴⁷³YSASSTA⁴⁸⁰
 Swiss-Prot-Link: P31749
 Akt2: L⁴⁶¹LELDQRTHFPQFS⁴⁷⁴YSASIRE⁴⁸¹
 Swiss-Prot-Link: P31751
 Akt3: D⁴⁶¹NERRPHFPQFS⁴⁷²YSASGRE⁴⁷⁹
 Swiss-Prot-Link: Q9Y243

Akt Proteins

Akt1 (human) (rec.) (GST-tag)

BML-SE416-0005 5 µg
 BML-SE416-0020 20 µg
 Produced in insect cells. Full length active enzyme. **APPLICATION:** Kinetic and functional studies, phosphorylation of target substrates, drug screening.

Akt2 (human), (rec.)

ADI-PPK-411-Z 5 µg
 Active recombinant human full length Akt2 protein expressed by baculovirus in Sf9 insect cells.

Akt2 (human), (rec.) (GST-tag)

BML-SE247-0020 20 µg
 Produced in a baculovirus expression system. The ΔPH domain of human Akt2 (PKBβ; protein kinase B β) (aa 119-481) is fused at the N-terminus to a GST-tag. **APPLICATION:** Useful for kinetic, substrate, inhibitor studies and drug screening.

[T³⁰⁸E, S⁴⁷³D]Akt2 (human) (rec.) (GST-tag)

BML-SE248-0020 20 µg
 Produced in insect cells. ΔPH domain of human Akt2 (PKBβ; protein kinase B β) (aa 119-481) fused to a GST-tag. Constitutively active due to S⁴⁷³D and T³⁰⁸E mutations. **APPLICATION:** Useful for kinetic, substrate and inhibitor studies. 5x more active than BML-SE247.

Akt3 (human), (rec.) (GST-tag)

BML-SE369-0005 5 µg
 BML-SE369-0020 20 µg
 Produced in insect cells. Full length active enzyme. **APPLICATION:** Kinetic and functional studies, phosphorylation of target substrates, drug screening.

Akt Substrates

Akt substrate

[H-Arg-Pro-Arg-Ala-Ala-Thr-Phe-OH]
 BML-P129-0001 1 mg
 Derived from variations on the GSK-3 phosphorylation site, it conforms to the Akt/PKB consensus motif, Arg-Xaa-Arg-Xaa-Xaa-(Ser/Thr). Unlike other Akt-PKB peptide substrates, it is not phosphorylated by p70 S6 kinase or MAPKAP kinase-1.

PAK4 & Akt substrate

[H-Cys-Lys-Arg-Pro-Arg-Ala-Ala-Ser-Phe-Ala-Glu-OH]
 BML-P194-0001 1 mg
 Suitable substrate for PAK4, Akt, and PAK7.

Crosstide

[H-Gly-Arg-Pro-Arg-Thr-Ser-Ser-Phe-Ala-Glu-Gly-OH]
 BML-P149-0001 1 mg
 This peptide sequence corresponds to the sequence of GSK-3 surrounding the serine phosphorylated by MAP kinase activated protein kinase (MAPKAP kinase-1) and p70 S6 kinase. Acts as a substrate for Akt (PKB), K_m=4µM. Useful in phosphocellulose kinase assays.

Crosstide (biotinylated)

[Biotin-Gly-Arg-Pro-Arg-Thr-Ser-Ser-Phe-Ala-Glu-Gly-OH]
 BML-P191-0001 1 mg
 This peptide sequence corresponds to the sequence of GSK-3 surrounding the serine phosphorylated by MAP kinase activated protein kinase (MAPKAP kinase-1) and p70 S6 kinase. Acts as a substrate for Akt (PKB), K_m=4µM. The N-terminal biotin allows this peptide to be used in kinase assays with streptavidin-bound membranes.

Akt Antibodies

Product	Specificity	Application	Prod. No.	Size
Akt, mAb (5C10)	Human, mouse, rat, dog	WB	ALX-804-553-C100	100 µg
Akt, mAb (19G7)	Human, mouse, rat, dog, hamster	ELISA, IP, WB	ALX-803-303-C100	100 µg
Akt, pAb	Human, mouse, rat, <i>drosophila</i>	WB	ALX-210-226-C200	200 µg
Akt, pAb	Human, mouse and rat	WB	ADI-KAP-PK004-D ADI-KAP-PK004-F	50 µg 200 µg
Akt (mouse), pAb	Mouse	ELISA, IP, WB	ALX-210-534-R100	100 µl
Akt (dephosphorylated), mAb (11A11)	Human, mouse, rat, dog	IHC, WB	ALX-804-400-C100	100 µg
[pSer⁴⁷³]Akt, mAb (11E6)	Human, mouse, rat, dog	FC, IHC, WB	ALX-804-401-C100	100 µg
[pSer⁴⁷³]Akt, pAb	Mouse, (human, rat, bovine and chicken)	WB	ADI-905-773-100	100 µg
[pThr³⁰⁸]Akt, pAb	Human and mouse	WB	ADI-KAP-PK007-E	100 µl
Akt1, mAb (5G12)	Human, rat, dog	IHC, WB	ALX-804-633-C100	100 µg
Akt1/2, pAb	Human, mouse and rat	WB	ADI-905-755-100	100 µg
Akt2, pAb	Human, mouse, rat, bovine, chicken, dog, guinea pig, hamster, monkey, pig, rabbit, sheep and <i>Xenopus</i>	ELISA, WB	ADI-KAP-PK008-E	100 µg
Akt2 (human), mAb (8B7)	Human	IHC, WB	ALX-804-634-C100	100 µg
Akt3 (human), mAb (66C1247.1)	Human	WB	ALX-804-379-C100	100 µg
Akt3, mAb (9B2)	Human and mouse	WB	ADI-905-821-100	100 µl

Akt Assay Kits

NEW

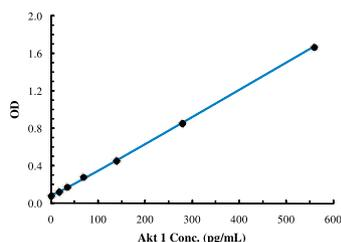
[pSer⁴⁷³]Akt1/2 EIA kit

ADI-900-162

96 wells

For the quantitative determination of [pSer⁴⁷³]Akt1/2 in cell lysates of human, mouse, and rat origin.

SENSITIVITY: 5.5 pg/ml (range 17.5 - 560 pg/ml).



Akt Kinase activity kit

ADI-EKS-400A

96 wells

For the measurement of Akt in partially purified, purified, or crude enzyme preparations from any species. This is a non-radioactive Akt kinase activity assay, providing a safe, rapid and reliable method for the screening of inhibitors or activators of Akt and for quantitating the activity of Akt in purified or partially purified enzyme preparations. The Akt Kinase activity kit is based on a solid phase enzyme immunoassay (EIA) that utilizes a specific synthetic peptide as a substrate for Akt and a polyclonal antibody that recognizes the phosphorylated form of the substrate. The assay is designed for the analysis of Akt (1, 2, 3) activity in the solution phase.

HIGHLIGHT

Akt Inhibitors

BML-257

BML-EI336-0010

10 mg

BML-EI336-0050

50 mg

Deguelin

ALX-350-118-M005

5 mg

ALX-350-118-M025

25 mg

(-)-Epigallocatechin gallate

ALX-270-263-M010

10 mg

ALX-270-263-M050

50 mg

ML-9

BML-EI153-0010

10 mg

BML-EI153-0050

50 mg

PKC-412

BML-EI370-0001

1 mg

BML-EI370-0005

5 mg

SH-5

ALX-270-349-MC05

0.5 mg

ALX-270-349-M001

1 mg

SH-6

ALX-270-350-MC05

0.5 mg

ALX-270-350-M001

1 mg

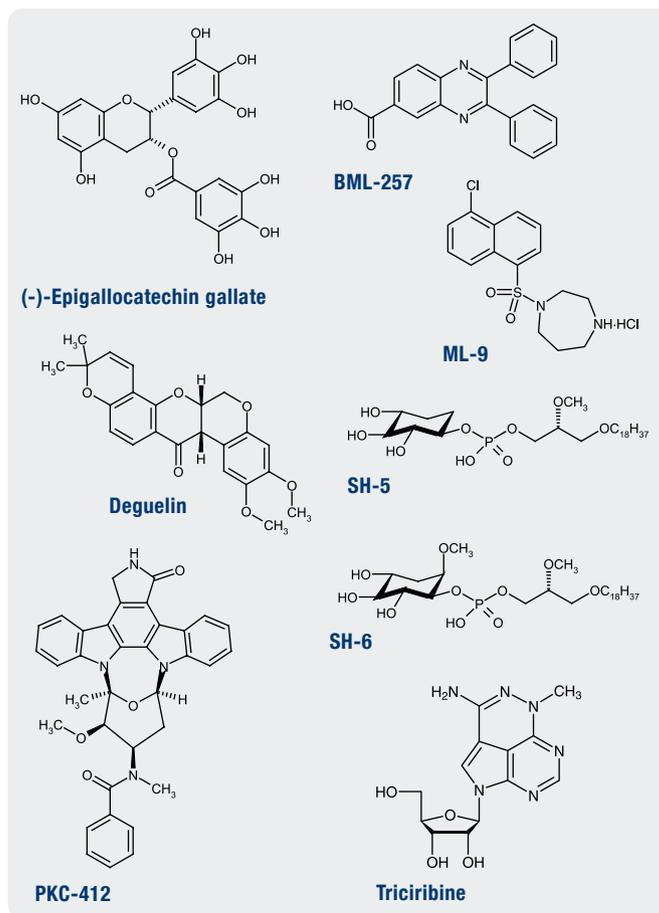
Triciribine

BML-EI332-0001

1 mg

BML-EI332-0005

5 mg



PDK

PDK1, pAb

ADI-KAP-PK112-D

50 µg

ADI-KAP-PK112-F

200 µg

From rabbit. **IMMUNOGEN:** Synthetic peptide derived from the sequence of human PDK1. **SPECIFICITY:** Recognizes human, mouse, rat, dog, monkey, pig and rabbit PDK1 (Pyruvate dehydrogenase kinase). **APPLICATION:** IHC, WB.

PDK1 (human), (rec.) (His-tag)

BML-SE351-0005

5 µg

BML-SE351-0020

20 µg

Produced in insect cells. Active, full length PDK1. **APPLICATION:** Kinetic and functional studies, phosphorylation of target substrates, drug screening.

PDKtide

BML-P250-0100

100 µg

This peptide, KTFCGTPEYLAPEVRREPRILSEEEQEMFRDFDYADWC, is derived from AKT1 and PKN2/PRK2. It is a substrate for PDK1, CDK9/Cyclin T1 and JAK2.

PDKtide (biotinylated)

BML-P251-0100

100 µg

This peptide, Biotin-Ahx-KTFCGTPEYLAPEVRREPRILSEEEQEMFRDFDYADWC, is derived from AKT1 and PKN2/PRK2. It is a substrate for PDK1, CDK9/Cyclin T1 and JAK2. The biotin allows peptide to be used in kinase assays with streptavidin-bound membranes.

SGK

SGK1 (human), (rec.) (GST-tag)

BML-SE398-0005	5 µg
BML-SE398-0020	20 µg

Produced in insect cells. **APPLICATION:** Kinetic and functional studies, phosphorylation of target substrates, drug screening.

SGK-1 (truncated NT 60 aa) (human), (rec.)

ADI-PPK-459-Z	5 µg
---------------	------

Active recombinant human SGK-1 containing an N-terminal GST tag, (truncated NT 60 a.a.).

SGK-1 (CT), pAb

ADI-KAP-PK015-C	25 µg
ADI-KAP-PK015-E	100 µg

From rabbit. **IMMUNOGEN:** Synthetic peptide derived from sequence near the C-terminus of human SGK-1 or OVA. **SPECIFICITY:** Recognizes human and monkey SGK-1. **APPLICATION:** IHC, WB.

SGK-1 (NT), pAb

ADI-KAP-PK016-C	25 µg
ADI-KAP-PK016-E	100 µg

From rabbit. **IMMUNOGEN:** Synthetic peptide derived from sequence near the N-terminus of human SGK-1 or OVA. **SPECIFICITY:** Recognizes human SGK-1. **APPLICATION:** WB.

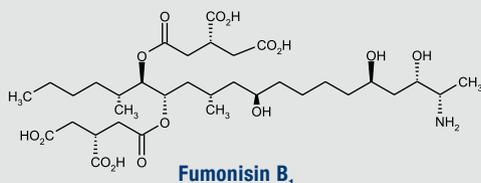
Akt Activator

Fumonisin B₁

BML-SL220-0001	1 mg
BML-SL220-0005	5 mg

A fungal metabolite produced by *Fusarium moniliforme*. Mycotoxin. Activates Akt (protein kinase B; PKB), which leads to increased survival, inhibition of GSK-3β activity and post-translational stabilization of cyclin D1. Inhibits sphinganine N-acyl-transferase (ceramide synthase) ($IC_{50}=0.1\mu\text{M}$). Inhibits de novo sphingolipid biosynthesis ($IC_{50}=0.7\mu\text{M}$ for sphingomyelin labeling by [³H]sphinganine, cultured cerebellar neurons). Inhibits growth of axons in cultured hippocampal neurons by blocking glycosphingolipid synthesis. Induces DNA damage in the liver and cancer. Induces apoptosis.

LIT: A potential mechanism for fumonisin B(1)-mediated hepatocarcinogenesis: cyclin D1 stabilization associated with activation of Akt and inhibition of GSK-3beta activity: D. Ramljak, et al.; *Carcinogenesis* **21**, 1537 (2000) • For a comprehensive bibliography please visit our website.



SGK2 (human), (rec.) (GST-tag)

BML-SE350-0005	5 µg
BML-SE350-0020	20 µg

Produced in insect cells. Active, full length serum- and glucocorticoid-induced kinase 2 is fused at the N-terminus to a GST-tag. **APPLICATION:** Kinetic and functional studies, phosphorylation of target substrates, drug screening.

SGK3 (human), (rec.) (GST-tag)

BML-SE415-0005	5 µg
BML-SE415-0020	20 µg

Produced in insect cells. Full length active enzyme. **APPLICATION:** Kinetic and functional studies, phosphorylation of target substrates, drug screening.

SGKtide, (substrate)

ADI-SPK-110-J	1 mg
---------------	------

Synthetic SGKtide substrate peptide.

PTEN

PTEN (human), (rec.) (His-tag)

BML-SE402-0010	10 µg
----------------	-------

Produced in *E. coli*. Full-length human PTEN (aa 1-403) fused to a N-terminal His tag.

PTEN, mAb (11G8.1)

ALX-804-451-C100	100 µg
------------------	--------

CLONE: 11G8.1. **ISOTYPE:** Mouse IgG. **IMMUNOGEN:** Recombinant human PTEN. **SPECIFICITY:** Recognizes human and mouse PTEN (epitope aa 152-200). Detects a band of ~60 kDa by Western blot. **APPLICATION:** IP, WB.

PTEN, mAb (6H2.1)

ALX-804-254-C100	100 µg
------------------	--------

CLONE: 6H2.1. **ISOTYPE:** Mouse IgG. **IMMUNOGEN:** Recombinant human PTEN. **SPECIFICITY:** Recognizes human, mouse and rat PTEN (epitope aa 304-403). Detects a band of ~60kDa by Western blot. **APPLICATION:** ELISA, IHC, ICC, IP, WB.

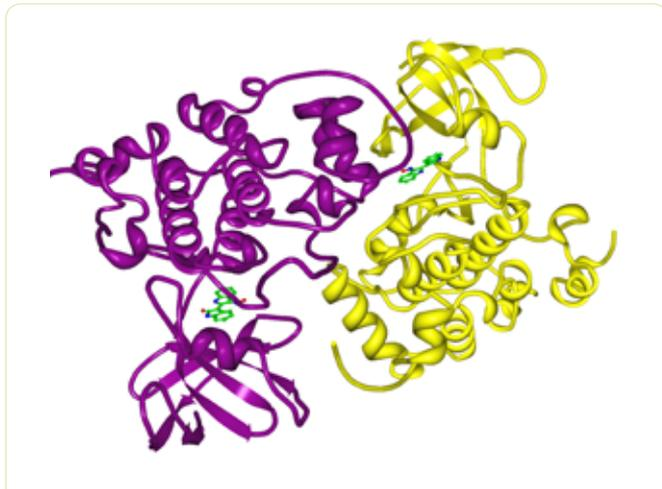
Ship2

Ship2 (catalytic domain) (human), (rec.)

BML-SE539-0010	10 µg
----------------	-------

Produced in *E. coli*. Recombinant catalytic domain of SHIP2 (Inpp1) fused to a His tag.

GSK-3 [Glycogen Synthase Kinase-3]



CRYSTAL STRUCTURE: Crystal structure of human GSK3-β complexed with Indirubin-3'-monoxime (Prod. No. ALX-270-271).

Glycogen synthase kinase-3 (GSK-3) is a multifunctional cytoplasmic serine/threonine kinase. This enzyme has several unique features compared to other kinases: i) GSK-3 is generally active ii) upon extracellular stimulation, GSK-3 gets inactivated through phosphorylation and iii) GSK-3 acts as a negative regulator in most pathways (except in NF-κB signaling).

GSK-3 is a key regulator of glycogen and protein synthesis and thus plays an important role in insulin signal transduction and metabolic regulation. Increased GSK-3 activity has been linked to pathology in diseases such as Alzheimer's disease and type 2 diabetes. Several GSK-3 inhibitors, such as the aloisines, the paullones and the maleimides, have been developed as therapeutic agents.

LITERATURE REFERENCES:

GSK-3: tricks of the trade for a multi-tasking kinase: B.W. Doble & J.R. Woodgett; *J. Cell Sci* **116**, 1175 (2003) (Review) • 9-cyano-1-azapallone (cazapallone), a glycogen synthase kinase-3 (GSK-3) inhibitor activating pancreatic beta cell protection and replication: H. Stukenbrock, et al.; *J. Med. Chem.* **51**, 2196 (2008) • Glycogen synthase kinase-3 (GSK-3) inhibitors reach the clinic: M. Medina & A. Castro; *Curr. Opin. Drug Discov. Devel.* **11**, 533 (2008) • Targeting glycogen synthase kinase-3 (GSK-3) in the treatment of Type 2 diabetes: K. MacAulay & J.R. Woodgett; *Expert Opin. Ther. Targets* **12**, 1265 (2008)

GSK-3 Protein

GSK-3β (human), (rec.) (GST-tag)

- BML-SE355-0005 5 µg
- BML-SE355-0020 20 µg

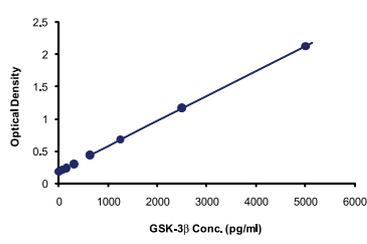
Produced in insect cells. Full length human GSK-3β is fused at the N-terminus to a GST-tag. **APPLICATION:** Kinetic and functional studies, phosphorylation of target substrates, drug screening.

GSK-3 ELISA Kits

NEW

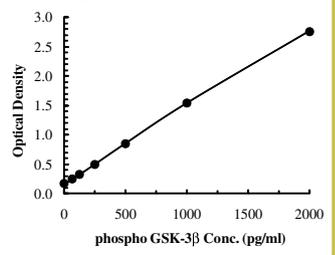
GSK-3β EIA kit

ADI-900-144 96 wells
For the quantitative determination of GSK-3β in cell lysates of human and rat origin. **SENSITIVITY:** 74.4 pg/ml (range 78.1 - 5,000 pg/ml).



[pSer⁹]GSK-3β EIA kit

ADI-900-123A 96 wells
For the quantitative determination of [pSer⁹]GSK-3β in cell lysates of human, mouse, and rat origin. **SENSITIVITY:** 9.0 pg/ml (range 62.5 - 2,000 pg/ml).



HIGHLIGHT

GSK-3 Antibodies

[pSer²¹]GSK-3 α , mAb (9B8)

ADI-905-759-100 100 μ g

CLONE: 9B8. **ISOTYPE:** Mouse IgG1. **IMMUNOGEN:** Synthetic phospho-peptide derived from the sequence of GSK-3 β , conjugated to KLH. **SPECIFICITY:** Recognizes human, mouse, rat and dog GSK-3 α . **APPLICATION:** ELISA, WB.

GSK-3 α / β , mAb (1H8)

ADI-KAM-ST002-D 50 μ g

ADI-KAM-ST002-F 200 μ g

CLONE: 1H8. **ISOTYPE:** Mouse IgG2b. **IMMUNOGEN:** Recombinant *Xenopus* GSK-3 β protein. **SPECIFICITY:** Recognizes human, mouse, rat, bovine, dog, hamster, monkey, pig, rabbit and sheep GSK-3 α / β . **APPLICATION:** WB.

GSK-3 α / β , mAb (1HB)

BML-SA364-0100 100 μ l

CLONE: 1HB. **ISOTYPE:** Mouse IgG2a. **IMMUNOGEN:** Recombinant *Xenopus laevis* glycogen synthase kinase-3 (GSK-3). **SPECIFICITY:** Recognizes human, mouse, rat and *X. laevis* GSK-3 α and β . **APPLICATION:** ELISA, IP, WB.

[pTyr²¹⁶/Tyr²⁷⁹]GSK-3 α / β , mAb (6D3)

ADI-905-762-100 100 μ g

CLONE: 6D3. **ISOTYPE:** Mouse IgG1. **IMMUNOGEN:** Synthetic phospho-peptide derived from the sequence of GSK, conjugated to hemocyanin. **SPECIFICITY:** Recognizes human and mouse GSK-3 α / β . **APPLICATION:** WB.

[pTyr²¹⁶/Tyr²⁷⁹]GSK-3 α / β , pAb

ADI-KAP-ST012-E 100 μ l

From rabbit. **IMMUNOGEN:** Human GSK-3 α / β synthetic peptide containing phosphorylated tyrosine 216/279. **SPECIFICITY:** Recognizes human and mouse GSK-3 α / β . **APPLICATION:** ELISA, WB.

GSK-3 β , mAb (11B9)

ALX-804-548-C100 100 μ g

CLONE: 11B9. **ISOTYPE:** Mouse IgG1. **IMMUNOGEN:** Synthetic peptide corresponding to the activation loop of GSK-3 β (glycogen synthase kinase-3 β). **SPECIFICITY:** Recognizes the activation loop of human, mouse, rat and dog GSK-3 β . **APPLICATION:** WB.

GSK-3 β , pAb

ADI-905-679-100 100 μ g

From rabbit. **IMMUNOGEN:** Synthetic peptide derived from sequence near the C-terminus of human GSK-3 β . **SPECIFICITY:** Recognizes human, mouse, rat and bovine GSK-3 β . **APPLICATION:** IHC, WB.

GSK-3 β , pAb

ADI-KAP-ST002-C 25 μ g

ADI-KAP-ST002-E 100 μ g

From rabbit. **IMMUNOGEN:** Synthetic peptide derived from the sequence of rat GSK-3 β ; sequence identical to human. **SPECIFICITY:** Recognizes human, mouse, rat and bovine GSK-3 β . **APPLICATION:** IHC, WB.

[pSer⁹]GSK-3 β , mAb (2D3)

ALX-804-374-C100 100 μ g

CLONE: 2D3. **ISOTYPE:** Mouse IgG1. **IMMUNOGEN:** Synthetic peptide corresponding to a portion of phosphorylated GSK-3 β (glycogen synthase kinase-3 β). **SPECIFICITY:** Recognizes human, mouse and dog GSK-3 β phosphorylated at Ser⁹. Does not cross-react with non-phosphorylated GSK-3 β , GSK-3 α or phosphorylated GSK-3 α . **APPLICATION:** WB.

[pSer⁹]GSK-3 β , mAb (3A8)

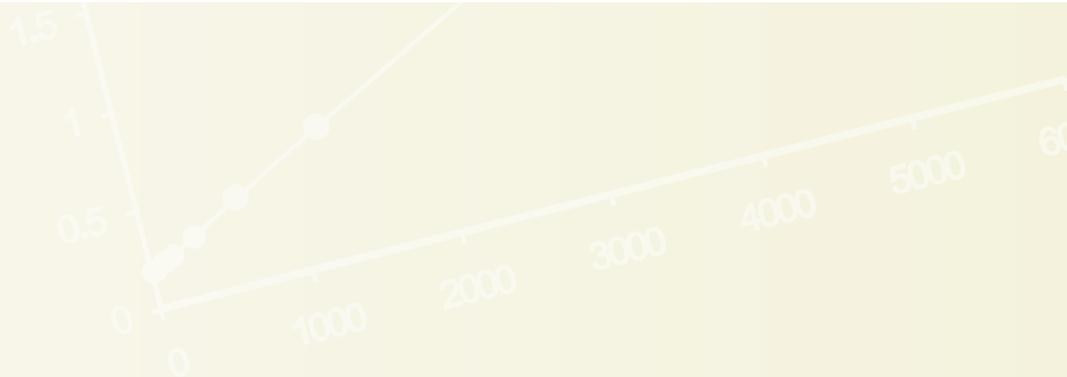
ADI-905-761-100 100 μ g

CLONE: 3A8. **ISOTYPE:** Mouse IgG1. **IMMUNOGEN:** Synthetic phospho-peptide derived from the sequence of GSK-3 β , conjugated to KLH. **SPECIFICITY:** Recognizes human, mouse, rat and dog GSK-3 β . **APPLICATION:** ELISA, WB.

[pSer⁹]GSK-3 β , pAb

ADI-KAP-ST011-E 100 μ g

From rabbit. **IMMUNOGEN:** Synthetic phospho-peptide derived from the sequence of human GSK-3 β . **SPECIFICITY:** Recognizes human, mouse, rat and *Xenopus* GSK-3 β . **APPLICATION:** ELISA, WB.



GSK-3 Inhibitors

Aloisine

[RP106]

ALX-270-386-M001 1 mg

Potent, cell permeable, selective ATP-competitive inhibitor of CDK1/cyclin B (IC_{50} =700nM), CDK5/p25 (IC_{50} =1.5 μ M), and GSK-3 (IC_{50} =920nM).

Aloisine A

[RP107]

ALX-270-385-M001 1 mg

Cell permeable, potent, selective, reversible and ATP-competitive inhibitor of CDK1/cyclin B (IC_{50} =150nM), CDK2/cyclin A (IC_{50} =120nM), CDK2/cyclin E (IC_{50} =400nM), CDK5/p25 (IC_{50} =200nM), CDK5/p35 (IC_{50} =160nM) and GSK-3 α (IC_{50} =500nM). Also inhibits GSK-3 β (IC_{50} =650nM) and c-Jun N-terminal kinase (JNK) (IC_{50} ~3-10 μ M).

Alsterpaullone

ALX-270-275-M001 1 mg

ALX-270-275-M005 5 mg

Potent inhibitor of CDK1/cyclin B (IC_{50} =35nM). Potent and selective inhibitor of CDK2/cyclin A, CDK2/cyclin E (IC_{50} =200nM), CDK5/p25 (IC_{50} =40nM), CDK5/p35 (IC_{50} =40nM) and GSK-3 β .

3-Amino-1H-pyrazolo[3,4-b]quinoxaline

ALX-270-387-M001 1 mg

ALX-270-387-M005 5 mg

Selective inhibitor of GSK-3 β (IC_{50} =1 μ M).

AR-A014418

ALX-270-468-M001 1 mg

Synthetic. Cell permeable, specific and potent inhibitor of GSK-3 (IC_{50} =104nM). Inhibition is competitive with respect to ATP (K_i =38nM).

1-Azakenpaullone

ALX-270-430-M001 1 mg

ALX-270-430-M005 5 mg

Potent and ATP-competitive inhibitor of GSK-3 β (IC_{50} =18nM).

6BIO

ALX-430-156-M001 1 mg

Potent, reversible and ATP-competitive inhibitor of GSK-3 α/β .

10Z-Hymenialdisine

ALX-350-289-C500 500 μ g

ALX-350-289-M001 1 mg

Isolated from sponge *Axinella carteri*. Inhibitor of DNA damage checkpoint at G2 phase (IC_{50} =6 μ M), cyclin-dependent kinases CDK1/cyclin B (IC_{50} =22nM), CDK2/cyclin A (IC_{50} =70nM), CDK2/cyclin E (IC_{50} =40nM), CDK4/cyclin D1 (IC_{50} =600nM), CDK5/p25 (IC_{50} =28nM), GSK-3 β (IC_{50} =10nM), and casein kinase 1 (CK1) (IC_{50} =35nM).

Indirubin

ALX-270-361-M001 1 mg

ALX-270-361-M005 5 mg

Synthetic. Inhibitor of cyclin-dependent kinases (CDK1/cyclin B (IC_{50} =10 μ M), CDK2/cyclin A (IC_{50} =2.2 μ M), CDK2/cyclin E (IC_{50} =7.5 μ M), CDK4/cyclin D1 (IC_{50} =12 μ M), CDK5/p35 (IC_{50} =5.5 μ M)) and of GSK-3 β (IC_{50} =600nM).

Indirubin-3'-monoxime

BML-CC207-0001 1 mg

BML-CC207-0005 5 mg

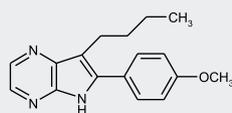
A potent inhibitor of glycogen synthase kinase-3 β (IC_{50} =22 nM). Also inhibits CDK1 (p34^{cdc2}; IC_{50} =180 nM) and CDK5 (IC_{50} =100 nM). It reversibly arrests asynchronous HBL-100 cells at G2. Induces apoptosis in the mammary carcinoma cell line MCF-7 (10 μ M).

Indirubin-5-sulfonic acid . sodium salt

ALX-270-296-M001 1 mg

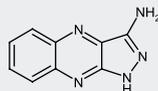
ALX-270-296-M005 5 mg

Inhibitor of cyclin-dependent kinases (CDKs), with selectivity for CDK1/cyclin B (IC_{50} =55nM), CDK2/cyclin A (IC_{50} =35nM), CDK2/cyclin E (IC_{50} =150nM), CDK4/cyclin D1 (IC_{50} =300nM), CDK5/p35 (IC_{50} =65nM) and GSK-3 β (IC_{50} =280nM).

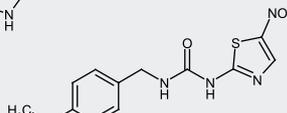
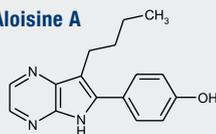


Aloisine

3-Amino-1H-pyrazolo[3,4-b]quinoxaline

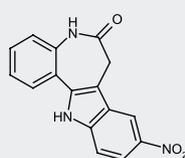
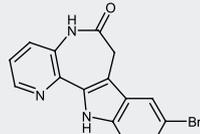


Aloisine A

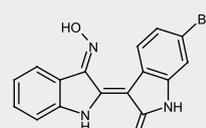


AR-A014418

1-Azakenpaullone

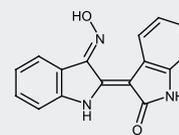
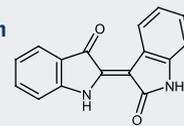


Alsterpaullone

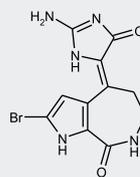


6BIO

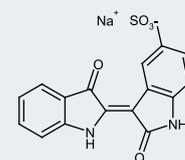
Indirubin



Indirubin-3'-monoxime



10Z-Hymenialdisine



Indirubin-5-sulfonic acid . sodium salt

5-Iodo-indirubin-3'-monoxime

ALX-270-424-M001 1 mg
 Potent inhibitor of GSK-3 β (IC_{50} =9nM).

Kenpaullone

BML-EI310-0001 1 mg
 BML-EI310-0005 5 mg
 Potent inhibitor of CDK1/cyclin B (IC_{50} =400nM). Also inhibits CDK2/cyclin A (IC_{50} =680nM), CDK5 (IC_{50} =850nM) and with much less effect other kinases. More recently, kenpaullone has been found to be a useful GSK-3 β inhibitor (IC_{50} =23nM).

Manzamine A

ALX-350-294-M001 1 mg
 Isolated from *Xestospongia sp.* GSK-3 Inhibitor.

SB-216763

BML-EI312-0001 1 mg
 BML-EI312-0005 5 mg
 Potent and selective inhibitor of GSK-3 β (IC_{50} =34nM). Stimulates glycogen synthesis in human liver cells and mimics other actions of insulin.

SB-415286

BML-EI311-0001 1 mg
 BML-EI311-0005 5 mg
 Potent and selective inhibitor of GSK-3 β (IC_{50} =78nM). Stimulates glycogen synthesis in human liver cells and mimics other actions of insulin.

Staurosporine

ALX-380-014-C100 100 μ g
 ALX-380-014-C250 250 μ g
 ALX-380-014-M001 1 mg
 ALX-380-014-M005 5 mg

Isolated from *Streptomyces staurosporeus*. Potent cell permeable inhibitor of a variety of protein kinases, e.g. GSK-3 β (IC_{50} =15nM).

Staurosporine

BML-EI156-0100 100 μ g
 BML-EI156-1000 1 mg

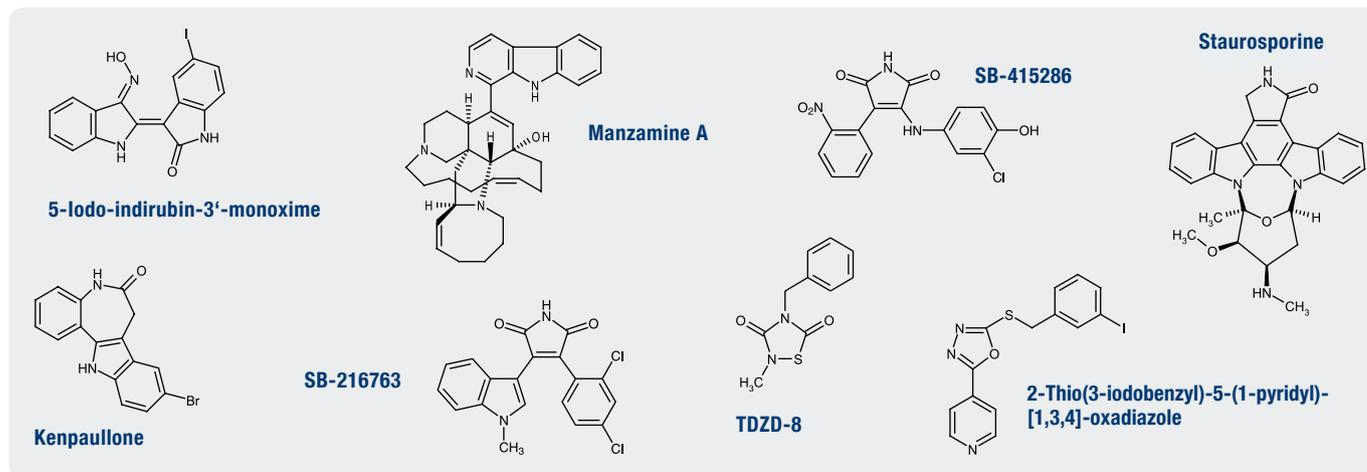
TDZD-8

ALX-270-354-M005 5 mg

Highly selective, non-ATP competitive inhibitor of GSK-3 β (IC_{50} =2 μ M). Binds to the active site of GSK-3 β . Does not significantly affect the activities of Cdk 1/cyclin B, CK-II, PKA, and PKC (IC_{50} >100 μ M).

2-Thio(3-iodobenzyl)-5-(1-pyridyl)-[1,3,4]-oxadiazole

ALX-270-358-M005 5 mg
 Potent inhibitor of GSK-3 β (IC_{50} =390nM).



GSK-3 Substrates

GSK-3 substrate

[H-Arg-Arg-Arg-Ala-Ala-Glu-Glu-Leu-Asp-Ser-Arg-Ala-Gly-pSer-Pro-Gln-Leu-OH]

BML-P151-0001 1 mg
 Corresponds to the sequence of the ϵ subunit of eIF2B surrounding serine 540, a target of GSK-3 (glycogen synthase kinase-3) phosphorylation, with the addition of a phosphoserine in the 4 position. This phosphoserine allows for more efficient phosphorylation of the second serine in the peptide (K_m =49 μ M for GSK-3 β , 83 μ M for GSK-3 α). Useful in phosphocellulose GSK-3 α and - β assays.

GSK-3 substrate II

[H-Tyr-Arg-Arg-Ala-Ala-Val-Pro-Pro-Ser-Pro-Ser-Leu-Ser-Pro-Ala-Ser-Ser-Pro-His-Gl n-pSer-Glu-Asp-Glu-Glu-OH]

BML-P193-0001 1 mg
 This peptide was derived from glycogen kinase I.

Nutrients & Nutrient Sensing

Nutrients (e.g. glucose, amino acids and fatty acids) can serve as important signaling molecules in complex pathways. Three major nutrient signaling pathways have been identified so far: the hexosamine signaling pathway, the mTOR signaling pathway and the AMPK pathway. These pathways regulate different aspects of energy metabolism and influence various cellular processes such as cell growth, proliferation and survival.

The hexosamine signaling pathway converts fructose-6-phosphate into uridine diphosphate N-acetyl glucosamine (UDP-GlcNAc). UDP-GlcNAc is a substrate for the addition (via an O-linkage) of a single N-acetylglucosamine to serine or threonine residues of nuclear and cytoplasmic proteins (O-glycosylation). O-glycosylation is a highly dynamic posttranslational modification that plays a role in key biological processes such as gene expression, cell growth, cell differentiation and fuel metabolism. Several studies have described a link between O-glycosylation and the induction of insulin resistance and type 2 diabetes. In addition, the hexosamine biosynthesis pathway regulates the secretion of both adiponectin and leptin from adipose tissue.

LITERATURE REFERENCE:
The hexosamine signaling pathway: deciphering the "O-GlcNAc code": D.C. Love & J.A. Hanover; *Sci. STKE* **2005**, re13

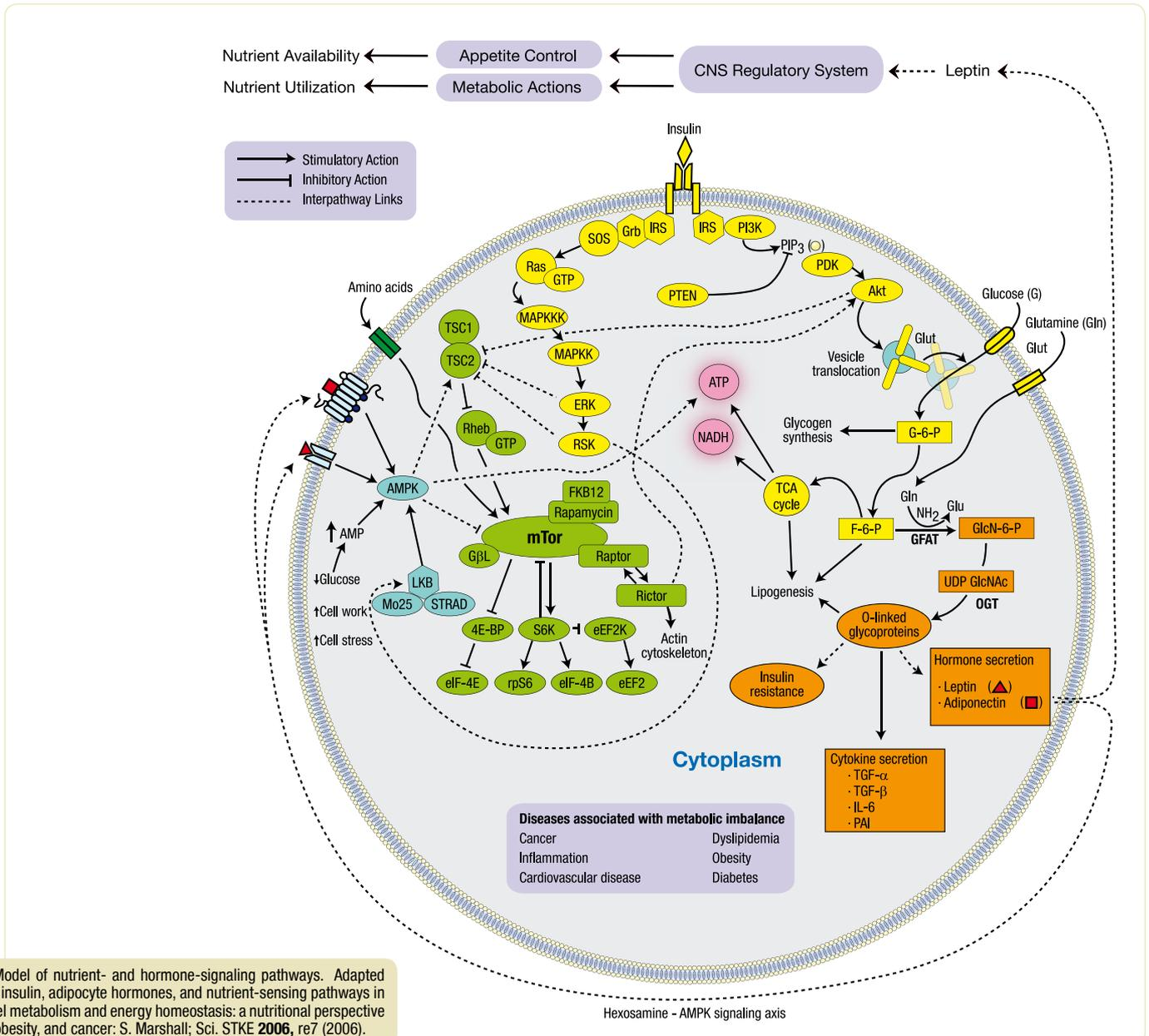


FIGURE 6: Model of nutrient- and hormone-signaling pathways. Adapted from: Role of insulin, adipocyte hormones, and nutrient-sensing pathways in regulating fuel metabolism and energy homeostasis: a nutritional perspective of diabetes, obesity, and cancer: S. Marshall; *Sci. STKE* **2006**, re7 (2006).

The mTOR Pathway

The mTOR (mammalian target of rapamycin) signaling pathway monitors intracellular amino acid availability and cellular energy status and links this information with external signals from cell surface receptors (such as the insulin receptor). This sensory input is then biochemically integrated and elicits a coordinated response that controls fundamental biological processes such as cell growth, cell proliferation and survival.

LITERATURE REFERENCE:

Expanding mTOR signaling: Q. Yang & K.L. Guan; Cell Res. **17**, 666 (2007)

mTOR (human FRB Domain), pAb

ALX-215-065-1

1 Vial

From rabbit. **IMMUNOGEN:** Recombinant human mTOR (FKBP12-rapamycin-binding (FRB) domain). **SPECIFICITY:** Recognizes the FRB domain of fusion proteins expressed in *Saccharomyces cerevisiae*. Not tested, but expected to recognize endogenous human FRB domain. **APPLICATION:** ICC, WB.

LT: The anchor-away technique: rapid, conditional establishment of yeast mutant phenotypes: H. Haruki, et al.; Mol. Cell **31**, 925 (2008)

mTOR, pAb

ALX-210-065-C050

50 µg

From rabbit. **IMMUNOGEN:** Human mTOR (mammalian target of rapamycin) (aa 1223-1290). **SPECIFICITY:** Recognizes human and rat mTOR. **APPLICATION:** IP, WB.

LT: Identification of TOR signaling complexes: more TORC for the cell growth engine: R.T. Abraham; Cell **111**, 9 (2002)

mTOR, pAb

ADI-905-687-100

100 µg

From rabbit. **IMMUNOGEN:** Synthetic peptide derived from the sequence of human mTOR. **SPECIFICITY:** Recognizes human and mouse mTOR. **APPLICATION:** WB.

Raptor, mAb (10E10)

ADI-905-765-100

100 µg

CLONE: 10E10. **ISOTYPE:** Mouse IgG1. **IMMUNOGEN:** Synthetic peptide derived from sequence of raptor, conjugated to hemocyanin. **SPECIFICITY:** Recognizes human, mouse, rat and dog raptor. **APPLICATION:** WB.

Rictor, mAb (1G11)

ADI-905-766-100

100 µg

CLONE: 1G11. **ISOTYPE:** Mouse IgG1. **IMMUNOGEN:** Synthetic phospho-peptide derived from the sequence of rictor, conjugated to hemocyanin. **SPECIFICITY:** Recognizes human, mouse and rat rictor. **APPLICATION:** WB.

Related Products

Rapamycin

BML-A275-0001

1 mg

BML-A275-0005

5 mg

Macrocyclic-triene antibiotic possessing potent immunosuppressant activity. It forms a complex with FKBP12 that binds to an effector, thus inhibiting IL-2 and other growth promoting lymphokines. The effectors were recently identified as FRAP (FKBP12 rapamycin-associated protein) and RAFT1 (rapamycin and FKBP12 target). Rapamycin/FKBP complex does not inhibit FRAP PI 4-kinase activity, but does inhibit FRAP autophosphorylation. Rapamycin induces inhibition of p70^{sgk}, p33^{cdk2} and p34^{cdc2}. Selectively blocks signaling leading to the activation of p70/85 S6 kinase. Enhances apoptosis. Activator of autophagy both *in vitro* and *in vivo*.

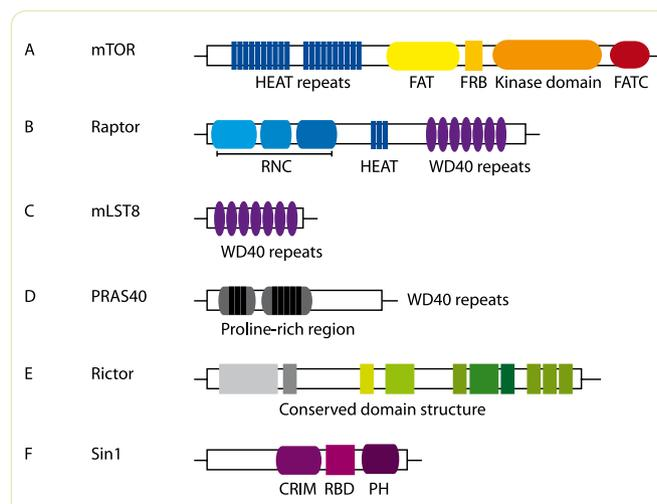


FIGURE 7: Schematic of mTOR complex components. Abbreviations: HEAT: a protein-protein interaction structure of two tandem anti-parallel ahelices found in huntingtin, elongation factor 3, PR65/A and TOR; FAT: α domain structure shared by FRAP, ATM and TRRAP; FRB: FKBP12/rapamycin binding domain; FATC: FAT C-terminus; RNC: Raptor N-terminal conserved domain; WD40: about 40 amino acids with conserved W and D forming four anti-parallel beta strands; CRIM: conserved region in the middle; RBD: Ras binding domain. Adapted from: Expanding mTOR signaling: Q. Yang & K.L. Guan; Cell Res. **17**, 666 (2007).

TSC1, pAb

ADI-905-713-100

100 µg

From rabbit. **IMMUNOGEN:** Synthetic peptide derived from the sequence of human TSC1. **SPECIFICITY:** Recognizes human, mouse and rat TSC1. **APPLICATION:** ICC, WB.

PRAS40 (human) (rec.)

BML-SE308-0025

25 µg

Produced in *E. coli*. Human PRAS40 (proline-rich Akt substrate 40) (aa 2-256).

[pThr²⁴⁶]PRAS40 (human), pAb

BML-SA360-0020

20 µl

BML-SA360-0100

100 µl

From rabbit. **IMMUNOGEN:** Synthetic peptide corresponding to a portion of human PRAS40 (proline-rich Akt substrate 40) phosphorylated at Thr²⁴⁶. **SPECIFICITY:** Recognizes PRAS40 phosphorylated at Thr²⁴⁶. **APPLICATION:** WB.

FKBP12, pAb

ALX-210-142-R100

100 µl

From rabbit. **IMMUNOGEN:** Synthetic peptide corresponding to aa 1-13 (G¹VQVETISP GDGR¹³) of human FKBP12 (FK506 binding protein 12kDa). **SPECIFICITY:** Recognizes human, mouse and rat FKBP12. **APPLICATION:** IHC (FS, PS), IP, WB. **BP:** ALX-156-002.

LT: The effects of FK506 on retinal ganglion cells after optic nerve crush: E.E. Freeman and C.L. Grosskreutz; IOVS **41**, 1111 (2000) • Immunophilin expression in the HIV-infected brain: C.L. Achim, et al.; J. Neuroimmunol. **157**, 126 (2004)

p70 S6K (substrate)

ADI-SPK-109-J 1 mg
Synthetic p70S6K/Rsk substrate peptide for use in kinase assays.

p70 S6K (human), (rec.)

BML-SE345-0005 5 µg
BML-SE345-0020 20 µg
Produced in insect cells. Full length active p70 S6 K (ribosomal protein S6 kinase (70kDa)).

[pThr³⁸⁹]p70 S6K (human), pAb

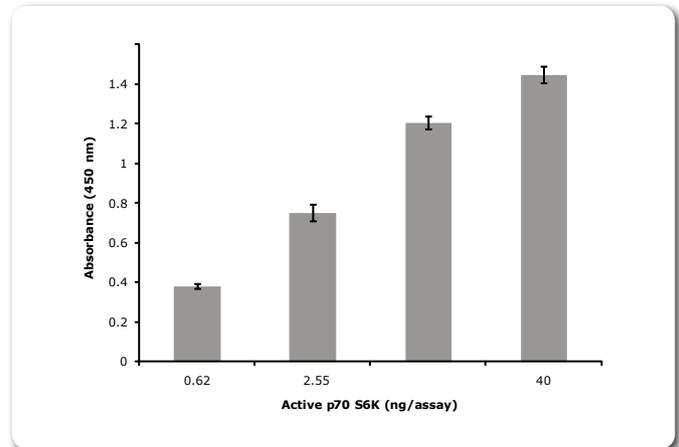
ALX-210-068-C050 50 µg
From rabbit. **IMMUNOGEN:** Synthetic peptide corresponding to aa 383-395 (CV³⁸³QFLGFpTYVAPSV³⁹⁵) of C-terminal human p70 S6K (ribosomal protein S6 kinase (70kDa)) phosphorylated at Thr³⁸⁹. **SPECIFICITY:** Recognizes human p70 S6K phosphorylated at Thr³⁸⁹. **APPLICATION:** WB.

p70 S6K, pAb

ADI-KAP-CC035-E 100 µg
From rabbit. **IMMUNOGEN:** Synthetic peptide derived from sequence near the C-terminus of human p70 S6K. **SPECIFICITY:** Recognizes human, mouse, rat and bovine p70 S6K. **APPLICATION:** ELISA, IP, WB.

p70 S6K activity kit

ADI-EKS-470 96 wells
For the measurement of p70 S6K activity in partially purified, purified, or crude enzyme preparations from any species.



The AMPK Pathway

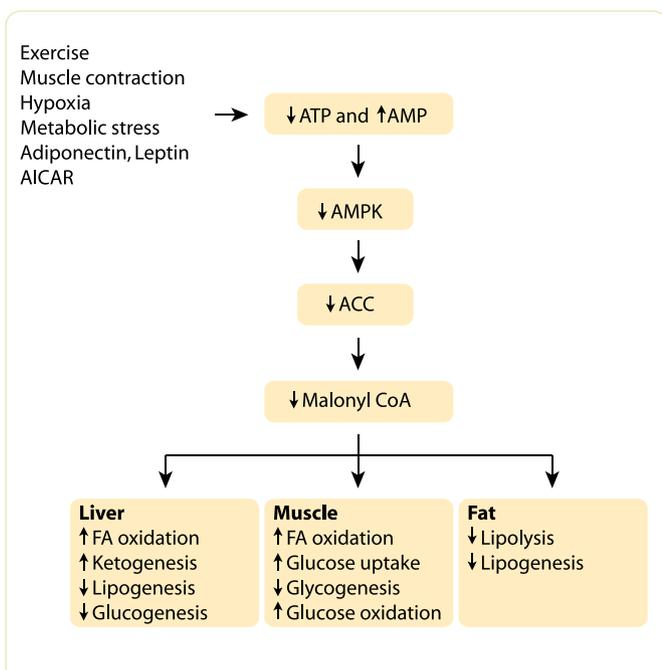


FIGURE 8: Metabolic effects of adenosine monophosphate activated protein kinase (AMPK). Abbreviations: AICAR = 5-aminoimidazole-4-carboxamide-ribofuranoside; ACC = Acetyl CoA carboxylase. Adapted from: Treating insulin resistance: future prospects: C.J. Bailey; Diab. Vasc. Dis. Res. 4, 20 (2007).

The serine-threonine kinase AMPK (AMP-activated protein kinase) is activated in response to low cellular energy and responds by suppressing cell growth and biosynthetic processes. One mechanism of how AMPK senses cellular energy levels involves the allosteric activation of the kinase activity of AMPK. Low cellular energy is characterized by reduced ATP and elevated AMP levels. In this case AMP directly activates AMPK and triggers a signal transduction cascade that regulates the activity of various downstream targets including transcription factors, enzymes and other regulatory proteins. A prominent downstream target of AMPK is mTOR. The phosphorylation of mTOR plays an important role in restoring ATP levels by inhibiting energy-consuming processes such as cell growth and protein synthesis.

In addition to allosteric activation, the activity of AMPK can also be regulated through phosphorylation. The serine-threonine kinase LKB1 as well as the hormones adiponectin and leptin can influence the phosphorylation state of AMPK.

AMPK is a key regulator of glucose and lipid metabolism. In the liver, activation of AMPK results in enhanced fatty acid oxidation and decreased production of glucose, cholesterol, and triglycerides. Pharmacological activation of AMPK in vivo improves blood glucose homeostasis, cholesterol concentrations and blood pressure in insulin-resistant rodents. This makes AMPK an attractive pharmacological target for the treatment of type 2 diabetes.

LITERATURE REFERENCES:

The role of AMP kinase in diabetes: P. Misra & R. Chakrabarti; Indian J. Med. Res. 125, 389 (2007) • AMP-activated protein kinase: Role in metabolism and therapeutic implications: G. Schimmack, et al.; Diabetes Obes. Metab. 8, 591 (2006)

AMPK Proteins

AMPK (human), (rec.) (His-tag)

BML-SE491-0005 5 µg

BML-SE491-0020 20 µg

Produced in insect cells. Full length active AMPK composed of subunits α 1, β 1 and γ 1. **APPLICATION:** Kinetic and functional studies, drug screening.

AMPK Substrates

SAMS peptide

[H-His-Met-Arg-Ser-Ala-Met-Ser-Gly-Leu-His-Leu-Val-Lys-Arg-Arg-OH]

BML-P231-0001 1 mg

This peptide is derived from rat acetyl-CoA carboxylase and is known to be a good substrate for AMPK.

SAMS peptide (biotinylated)

[Biotin-His-Met-Arg-Ser-Ala-Met-Ser-Gly-Leu-His-Leu-Val-Lys-Arg-Arg-OH]

BML-P232-0001 1 mg

This peptide contains a biotin group at its N- terminus. It is derived from rat acetyl-CoA carboxylase and is known to be a good substrate for AMPK. The biotin allows peptide to be used in kinase assays with streptavidin-bound membranes.

SAMS peptide (phosphorylated) (biotinylated)

[Biotin-His-Met-Arg-Ser-Ala-Met-pSer-Gly-Leu-His-Leu-Val-Lys-Arg-Arg-OH]

BML-P233-0001 1 mg

This peptide contains a biotin group at its N- terminus and a phospho-serine. It is derived from rat acetyl-CoA carboxylase. It may be used as a control peptide for unphosphorylated SAMS Peptide (BML-P231 and BML-P232). The biotin allows peptide to be used in kinase assays with streptavidin-bound membranes.

AMARA peptide substrate

[H-Ala-Met-Ala-Arg-Ala-Ala-Ser-Ala-Ala-Ala-Leu-Ala-Arg-Arg-Arg-OH]

BML-P270-0001 1 mg

This peptide is a substrate for AMPK and SIK kinases.

AMPK Inhibitor

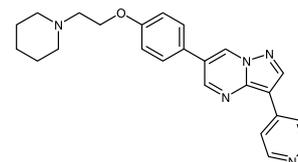
BML-275

[Compound C; Dorsomorphin]

BML-EI369-0005 5 mg

BML-EI369-0025 25 mg

Cell permeable pyrazolopyrimidine derivative that inhibits AMP kinase ($K_i=109\text{nM}$ in the absence of AMP) in an ATP-competitive manner. It displays no significant inhibition of ZAPK, SYK, PKCT, PKA and JAK3. Decreases food intake in mice and inhibits the effects of AICAR and metformin. It has recently been shown to inhibit BMP type I receptors ALK2, ALK3 and ALK6.



AMPK Activator

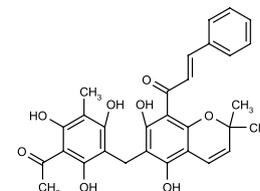
Rottlerin

[Mallotoxin]

ALX-350-075-M010 10 mg

ALX-350-075-M025 25 mg

Isolated from *Mallotus philippinensis*. Mitochondrial uncoupler that depolarizes the mitochondrial membrane potential, reduces cellular ATP levels, activates 5'-AMP-activated protein kinase (AMPK) and affects mitochondrial production of reactive oxygen species (ROS). Potent activator of multiple Ca^{2+} -sensitive K^+ channels. **lit:** The mouse ear edema: a quantitatively evaluable assay for tumor promoting compounds and for inhibitors of tumor promotion: M. Gschwendt, et al.; Cancer Lett. **25**, 177 (1984) • Rottlerin, a novel protein kinase inhibitor: M. Gschwendt, et al.; BBRC **199**, 93 (1994)



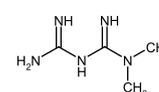
Metformin

ALX-270-432-G001 1 g

ALX-270-432-G005 5 g

Antidiabetic agent. Reduces the cardiovascular complications of diabetes. Does not induce hypoglycemia.

lit: Effects of metformin and thiazolidinediones on suppression of hepatic glucose production and stimulation of glucose uptake in type 2 diabetes: a systematic review: A. Natali and E. Ferrannini; Diabetologia **49**, 434 (2006) • For a comprehensive bibliography please visit our website.



Latest Insight

AMPK is strongly stimulated by exercise and plays a central role in sensing energy status. A recent study found that treatment of sedentary mice with AICAR mimicked many of the positive effects of exercise. Four weeks of AICAR treatment induced metabolic genes, decreased fat mass, increased oxygen consumption and enhanced running endurance by 44%. This study not only helps to characterize the molecular pathways that lead to the beneficial effects of exercise, but also suggests that exercise mimetic drugs may have therapeutic potential in treating muscle diseases like wasting and frailty or more broadly for treatment of obesity or heart disease where exercise has positive benefits.

lit: AMPK and PPARdelta agonists are exercise mimetics: V.A. Narkar et al.; Cell **134**, 405 (2008)

AICAR

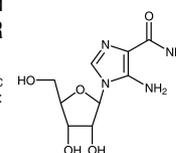
[5-Aminoimidazole-4-carboxamide 1- β -D-ribofuranoside]

BML-EI330-0050 50 mg

BML-EI330-0250 250 mg

Activates AMP Kinase in whole cells. Mimics the effects of insulin on the expression of two gluconeogenic genes PEPCK and glucose-6-phosphatase. Inhibits PPAR α coactivation and adipocyte differentiation. Substrate for the AICAR transformylase activity of ATIC.

lit: 5-aminoimidazole-4-carboxamide ribonucleoside. A specific method for activating AMP-activated protein kinase in intact cells?: J.M. Corton et al.; Eur. J. Biochem. **229**, 558 (1995)



PPARs [Peroxisome Proliferator-activated Receptor]

The PPAR (Peroxisome proliferator-activated receptor) family members play important roles in glucose and lipid metabolism and atherosclerosis. PPARs are ligand-activated transcription factors belonging to the superfamily of nuclear receptors. Both endogenous and pharmacologic ligands exist for PPARs. Upon binding, the ligands can heterodimerize with the retinoid X receptor, bind to the PPAR response element of its target gene and either activate or repress gene expression. Three forms of PPARs have been identified α , γ and β/δ . The three subtypes exist in different tissues in varying quantities with different effects (see Table below).

PPAR γ is highly expressed in adipose tissue and is involved in fatty acid uptake, glucose homeostasis and insulin resistance. It is the primary target of the thiazolidinediones (TZDs). The TZDs constitute a class of oral anti-diabetic agents that exert insulin-sensitizing effects in adipose tissue, skeletal muscle, and liver. In addition, TZDs also have beneficial effects on other components of the metabolic syndrome. They may influence cardiovascular risk and decrease inflammatory responses in insulin resistance. Within adipocytes, TZDs suppress the synthesis of IL-6, TNF- α , PAI-1, MCP-1 and angiotensinogen. Thus, TZDs have been used in patients with type 2 diabetes and pre-diabetic insulin resistance.

LITERATURE REFERENCES:

Peroxisome proliferator-activated receptor (PPAR) in metabolic syndrome and type 2 diabetes mellitus: M.A. Jay & J. Ren; *Curr. Diabetes Rev.* **3**, 33 (2007) (Review) • Inflammation in diabetes mellitus: role of peroxisome proliferator-activated receptor-alpha and peroxisome proliferator-activated receptor-gamma agonists: P. Libby & J. Plutzky; *Am. J. Cardiol.* **99**, 27B (2007) (Review) • Effects of peroxisome proliferator-activated receptors on lipoprotein metabolism and glucose control in type 2 diabetes mellitus: R.S. Rosenson; *Am. J. Cardiol.* **99**, 96B (2007) (Review) • Fat and beyond: the diverse biology of PPARgamma: P. Tontonoz & B.M. Spiegelman; *Annu. Rev. Biochem.* **77**, 289 (2008) (Review)

Subtype	Distribution	Physiological involvement
PPAR α	Liver, kidneys, heart, gut, skeletal muscle, adipose tissue	Lipid catabolism and oxidation, gluconeogenesis
PPAR β/δ	Ubiquitous	Adipocyte differentiation (minor), function incompletely understood
PPAR γ	Adipose tissue, large intestine, hematopoietic cells, kidney, liver, intestinal mucosa	Glucose and fatty acid uptake, gluconeogenesis, lipogenesis, glycogenesis, adipocytes differentiation, macrophage maturation, modulation of inflammation

TABLE: Distribution and actions of the PPARs. Adapted from: PPAR α : therapeutic role in diabetes-related cardiovascular disease: A.Y. Cheng & L.A. Leiter; *Diabetes Obes. Metab.* **10**, 691 (2008).

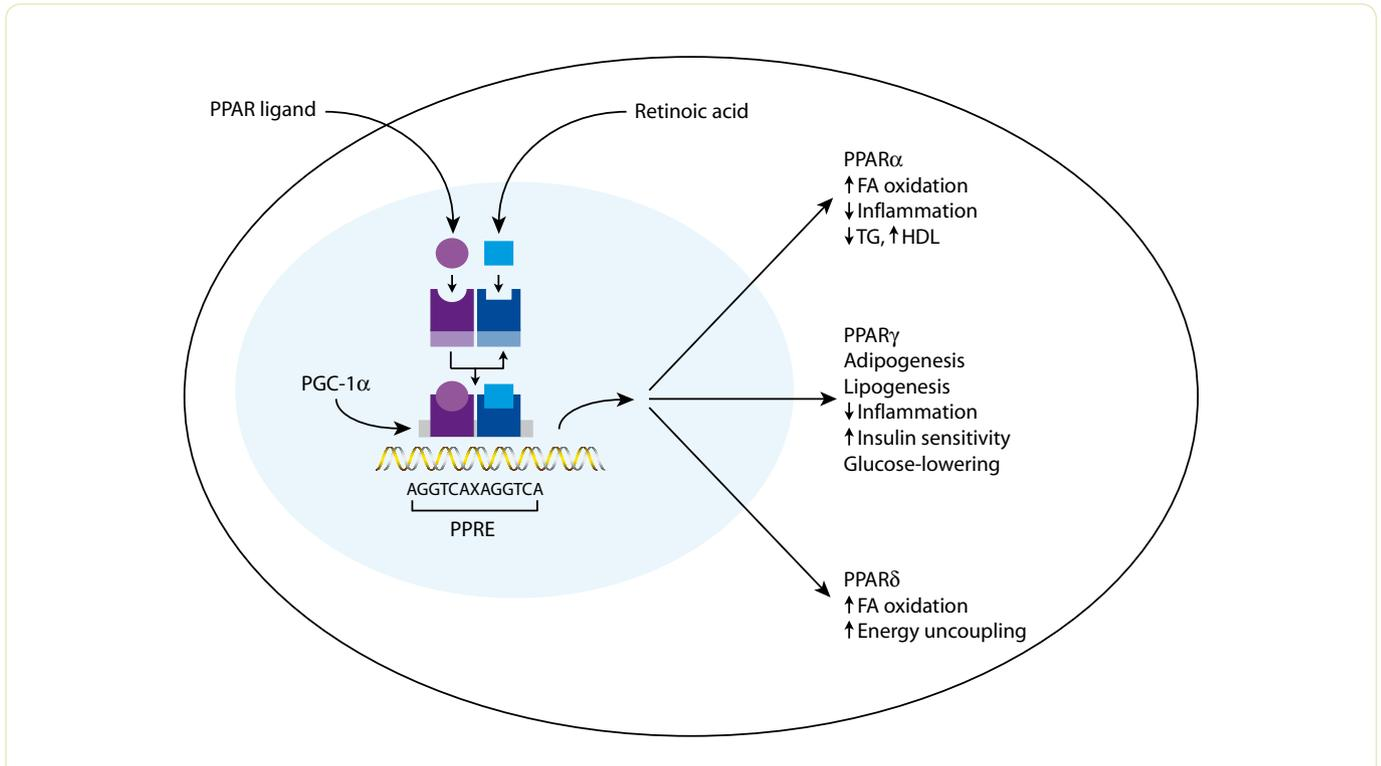


FIGURE 9: Actions of peroxisome proliferator-activated receptor (PPAR) ligands. Abbreviations: PGC-1 α = PPAR coactivator PGC-1 α ; FA = fatty acid; TG = triglyceride; HDL = high-density lipoprotein; PPRE = PPAR response element. Adapted from: *Treating insulin resistance: future prospects*: C.J. Bailey; *Diab. Vasc. Dis. Res.* **4**, 20 (2007).

PPAR Antibodies

PPAR, pAb

ALX-210-117-R100 100 µl

From rabbit. **IMMUNOGEN:** Synthetic peptide corresponding to aa 484-498 (¹⁴⁸⁴KKTETDMSLHPLLQ⁴⁹⁸) of the C-terminal domain of rat PPAR_γ2 (peroxisome proliferator activated receptor _γ2). This sequence is completely conserved in human, mouse, rat, bovine, chicken and pig PPAR_γ1, PPAR_α and NUC1. **SPECIFICITY:** Recognizes mouse and rat PPAR_γ2. Detects a band of ~55 kDa by Western blot. Cross-reacts with recombinant rat PPAR_α in Western blot and recombinant mouse PPAR_γ1 in immunoprecipitation. **APPLICATION:** IP, WB.

lit: mPPAR gamma 2: tissue-specific regulator of an adipocyte enhancer: P. Tontonoz, et al.; *Genes Dev.* **8**, 1224 (1994) • Differential expression of peroxisome proliferator-activated receptors (PPARs): tissue distribution of PPAR-alpha, -beta, and -gamma in the adult rat: O. Braissant, et al.; *Endocrinology* **137**, 354 (1996)

PPAR_α, mAb (3B6/PPAR)

ALX-804-255-C100 100 µg

CLONE: 3B6/PPAR. **ISOTYPE:** Mouse IgG2b. **IMMUNOGEN:** Purified recombinant PPAR_α (peroxisome proliferator activated receptor _α). **SPECIFICITY:** Recognizes human, mouse and rat PPAR_α. Detects a band of ~52 kDa by Western blot. **APPLICATION:** IP, WB.

lit: Evidence that PPARalpha is complexed with the 90 kDa heat shock protein and the hepatitis virus B X-associated protein 2: W.K. Sumanasekera, et al.; *J. Biol. Chem.* **278**, 4467 (2003) • PPARalpha activators inhibit vascular endothelial growth factor receptor-2 expression by repressing Sp-1-dependent DNA binding and transactivation: M. Meissner, et al.; *Circ. Res.* **94**, 324 (2004)

PPAR_α (mouse), pAb

ALX-210-190/1-R100 100 µl

From rabbit. **IMMUNOGEN:** Synthetic peptide corresponding to aa 1-18 (M¹VDTESPICPLSLEADD¹⁸) of mouse PPAR_α (peroxisome proliferator activated receptor _α). A single aa substitution (I8 to L8) exists between the mouse and human protein. **SPECIFICITY:** Recognizes mouse PPAR_α. Detects a band of ~52kDa by Western blot. Does not cross-react with PPAR_δ or PPAR_γ. **APPLICATION:** ICC, IP, WB. **BP:** ALX-165-029.

[pSer¹²]PPAR_α (mouse), pAb

ALX-210-361-C100 100 µg

From rabbit. **IMMUNOGEN:** Synthetic peptide corresponding to aa 8-19 (¹⁸CPLpSLEADDL¹⁹) of mouse PPAR_α (peroxisome proliferator activated receptor _α) phosphorylated at Ser¹². **SPECIFICITY:** Recognizes mouse PPAR_α phosphorylated at Ser¹². Detects a band of ~52 kDa by Western blot. **APPLICATION:** WB. **BP:** ALX-165-055.

lit: MCF-7 and T47D human breast cancer cells contain a functional peroxisomal response: M.W. Kilgore, et al.; *Mol. Cell Endocrinol.* **129**, 229 (1997) • Diverse signaling pathways modulate nuclear receptor recruitment of N-CoR and SMRT complexes: R.M. Lavinsky, et al.; *PNAS* **95**, 2920 (1998)

[pSer²¹]PPAR_α (mouse), pAb

ALX-210-362-C100 100 µg

From rabbit. **IMMUNOGEN:** Synthetic peptide corresponding to aa 14-25 (L¹⁴EADDLepSPLSE²⁵) of mouse PPAR_α (peroxisome proliferator activated receptor _α) phosphorylated at Ser²¹. **SPECIFICITY:** Recognizes mouse PPAR_α phosphorylated at Ser²¹. Detects a band of ~52kDa by Western blot. **APPLICATION:** WB. **BP:** ALX-165-056.

lit: See Prod. No. ALX-210-361.

PPAR_γ2 (mouse), pAb

ALX-210-192-R100 100 µl

From rabbit. **IMMUNOGEN:** Synthetic peptide corresponding to aa 1-16 (M¹GETLGDSPIDPESDS¹⁶) of human PPAR_γ2 (peroxisome proliferator activated receptor _γ2). **SPECIFICITY:** Recognizes mouse PPAR_γ2. Detects a band of ~56kDa by Western blot. Does not cross-react with PPAR_α or PPAR_δ. **APPLICATION:** ICC, WB. **BP:** ALX-165-031.

lit: Diverse signaling pathways modulate nuclear receptor recruitment of N-CoR and SMRT complexes: R.M. Lavinsky, et al.; *PNAS* **95**, 2920 (1998) • Delayed activation of PPARgamma by LPS and IFN-gamma attenuates the oxidative burst in macrophages: A.A. Von Knethen & B. Brune; *FASEB J.* **15**, 535 (2001) • Activation of peroxisome proliferator-activated receptor gamma by nitric oxide in monocytes/macrophages down-regulates p47phox and attenuates the respiratory burst: A. Von Knethen & B. Brune; *J. Immunol.* **169**, 2619 (2002) • A human peripheral blood monocyte-derived subset acts as pluripotent stem cells: Y. Zhao, et al.; *PNAS* **100**, 2426 (2003)

PPAR_δ, pAb

ALX-210-191/1-R100 100 µl

From rabbit. **IMMUNOGEN:** Synthetic peptide corresponding to aa 1-14 (M¹EQQEETPEAREE¹⁴) of mouse PPAR_δ (peroxisome proliferator activated receptor _δ). This sequence is ~86% conserved in human PPAR_δ. **SPECIFICITY:** Recognizes mouse and rat PPAR_δ. Detects a band of ~49kDa by Western blot. Does not cross-react with PPAR_α or PPAR_γ. **APPLICATION:** IHC (FS), ICC, WB. **BP:** ALX-165-030.



PPAR Modulators

BADGE

ALX-270-353-G025 25 g
Antagonist of PPAR γ . Binds to PPAR γ with an apparent Kd of 100 μ M.

Bezafibrate

BML-GR211-0001 1 g
BML-GR211-0005 5 g
Synthetic PPAR agonist. Induces fatty acid peroxisomal β -oxidation in rat white adipose tissue. Displays anti-inflammatory, anti-atherosclerosis and anti-obesity activity.

Carnosic acid

ALX-270-264-M010 10 mg
ALX-270-264-M050 50 mg
Isolated from *Rosmarinus officinalis*. Inhibits lipid peroxidation induced by NADH or NADPH oxidation. PPAR γ activator.

Ciglitazone

BML-GR205-0005 5 mg
BML-GR205-0025 25 mg
Synthetic. Selective PPAR γ agonist (EC_{50} =3 μ M).

Ciprofibrate

ALX-270-475-M025 25 mg
ALX-270-475-M100 100 mg
ALX-270-475-M250 250 mg
Hypolipidemic compound. PPAR α agonist.

Clofibrate

ALX-270-256-M500 500 mg
ALX-270-256-G001 1 g
Activates PPAR α and induces cytochrome P450 4A1 and 4A3.

5,8,11,14-Eicosatetraynoic acid

BML-ET004-0020 20 mg
BML-ET004-0100 100 mg

Inhibits arachidonic acid uptake. Inhibits cyclooxygenase (ID_{50} =8 μ M) and all lipoxygenases (ID_{50} =10, 0.3, and 0.2 μ M for 5-, 12- and 15-LO respectively) in whole cells. Inhibits PLA $_2$ and cytochrome P-450. Modulates Ca $^{2+}$ entry into cells. Stimulates luteinizing hormone release from cultured pituitary cells. PPAR agonist.

Fenofibrate

ALX-270-481-G005 5 g
Hypolipidemic compound. PPAR α agonist (EC_{50} =18 μ M for mouse and EC_{50} =30 μ M for human PPAR α). Also binds to PPAR γ , but with at least 10-fold less affinity and is inactive at PPAR δ (up to EC_{50} =100 μ M).

GW1929

ALX-420-029-M001 1 mg
ALX-420-029-M005 5 mg
Potent (<10nM) and subtype-selective (>1'000-fold) PPAR γ agonist which does not contain a thiazolidinedione moiety. The glucose-lowering effect of GW1929 in rats is 100-fold more potent than that of troglitazone (Prod. No. ALX-270-355).

GW501516

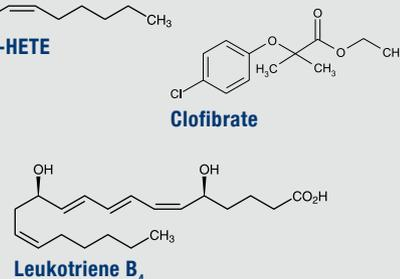
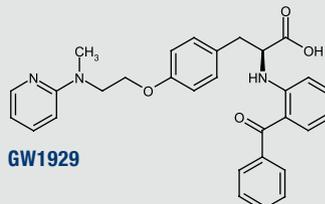
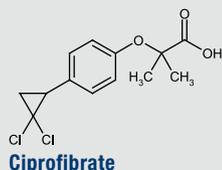
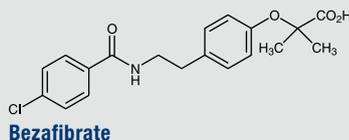
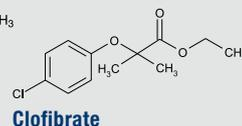
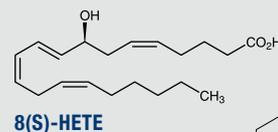
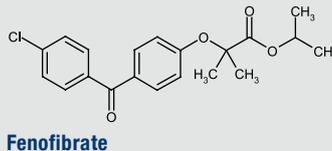
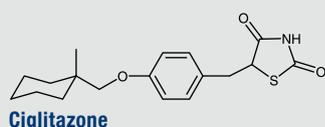
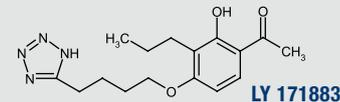
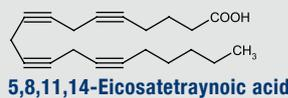
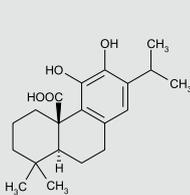
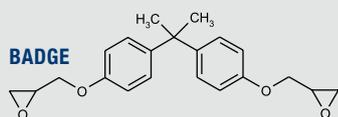
ALX-420-032-M001 1 mg
ALX-420-032-M005 5 mg
Specific agonist for PPAR δ (EC_{50} =1.1nM) with a 1'000-fold selectivity over the other human subtypes.

GW-9662

BML-GR234-0010 10 mg
BML-GR234-0050 50 mg
Blocks the PPAR γ -induced differentiation of monocytes to osteoclasts by >90% at a dose of 0.1 μ M. Much more potent antagonist than BADGE (Prod. No. ALX-270-353).

8(S)-HETE

BML-H008-0050 50 μ g
Activator of PPAR α .



Leukotriene B₄

BML-LB004-0050 50 µg
 Activator of PPAR α . Stimulates a number of leukocyte functions, including aggregation, stimulation of ion fluxes, enhancement of lysosomal enzyme release, superoxide anion production, chemotaxis and chemokinesis.

LY 171883

BML-RA101-0010 10 mg
 BML-RA101-0050 50 mg
 Competitive leukotriene D₄ (LTD₄) antagonist. Has also been shown to bind to the γ -isoform of the peroxisome proliferator activated receptor (PPAR γ).

MK-886

BML-EI266-0005 5 mg
 BML-EI266-0025 25 mg
 Inhibitor of PPAR α activation. Inhibits 5-lipoxygenase activating protein (FLAP) and leukotriene biosynthesis (IC₅₀=102nM).

Naproxen

ALX-270-102-G005 5 g
 PPAR agonist. Widely used non-steroidal anti-inflammatory drug.

N-Oleylethanolamide

BML-SL231-0010 10 mg
 BML-SL231-0050 50 mg
 Activates TRPV1. Does not activate cannabinoid receptors (CB) but is a PPAR α agonist (EC₅₀=120nM) *in vitro* and *in vivo*; induces satiety through activation of PPAR α . Inhibits ceramidase.

Pioglitazone

ALX-270-367-M001 1 mg
 ALX-270-367-M005 5 mg
 Pioglitazone selectively activates PPAR γ . It is about one tenth as potent as rosiglitazone (EC₅₀~500nM for human and mouse PPAR γ).

15-Deoxy- $\Delta^{12,14}$ -prostaglandin J₂

BML-PG050-0001 1 mg
 BML-PG050-0005 5 mg
 Suggested natural ligand for PPAR γ .

Pseudolaric acid B

ALX-350-108-MC01 0.1 mg
 ALX-350-108-M001 1 mg
 Isolated from *Pseudolarix kaempferi*. Activates PPAR α and the phospholipase C (PLC) signaling pathway. Stimulates peroxisomal fatty acyl-CoA oxidase activity.

Rosiglitazone

ALX-350-125-M025 25 mg
 ALX-350-125-M100 100 mg
 Potent insulin sensitizing agent. PPAR γ agonist. Hepatotoxic. Antidiabetic.

T0070907

BML-GR212-0010 10 mg
 BML-GR212-0050 50 mg
 Potent and selective PPAR γ antagonist. Blocks PPAR γ function in both cell-based reporter gene and adipocyte differentiation assays.

Tetradecylthioacetic acid

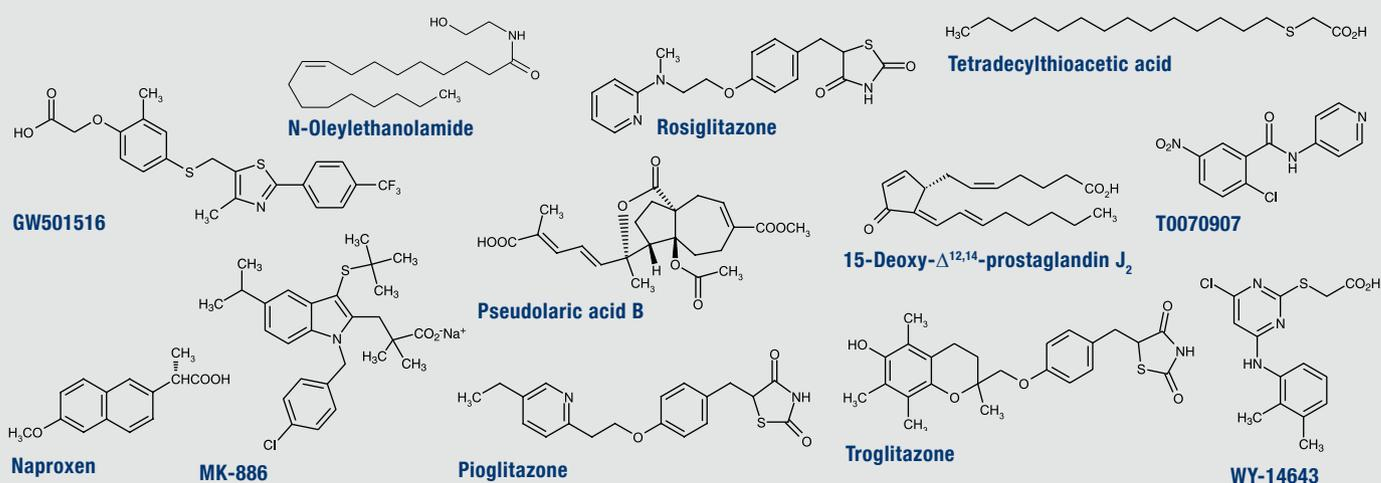
BML-GR236-0010 10 mg
 BML-GR236-0050 50 mg
 Tetradecylthioacetic acid (TTA) is a novel thio-fatty acid which cannot undergo β -oxidation. Activates rodent PPARs with rank order PPAR α > PPAR δ > PPAR γ .

Troglitazone

BML-GR210-0005 5 mg
 BML-GR210-0025 25 mg
 Synthetic. Potent and selective PPAR γ ligand. Binds to the PPAR γ ligand-binding domain and recruits ligand-specific transcriptional coactivators at 1µM. In skeletal muscle and adipocyte cell cultures, the EC₅₀ of troglitazone is about 10µM.

WY-14643

BML-GR200-0050 50 mg
 BML-GR200-0250 250 mg
 Widely used activator of PPARs.



Proteolytic Regulation of GLP-1 by DPPIV and Neprilysin

Glucagon like peptide-1 (GLP-1), along with glucose-dependent insulinotropic polypeptide (GIP), are incretin peptide hormones, which are released upon ingestion of food. Proglucagon is secreted and processed by proprotein convertases PC1/3 and PC2 into glucagon, GLP-1, GLP-2, and other bioactive peptides, depending on tissue type. GLP-1 (7-36) amide stimulates insulin and somatostatin secretion, and inhibits secretion of glucagon, ultimately resulting in satiety [1]. It is cleaved by dipeptidyl peptidase 4 (DPPIV) and neprilysin (NEP) into GLP-1 (9-36) amide, a form, that reverses the glucoregulatory actions of GLP-1 (7-36) amide [1-3]. Thus incretin mimetics, DPPIV/NEP inhibitors, and GLP-1 (9-36) amide peptides are of interest for treatment of diabetes and obesity [4]. Several DPPIV inhibitors are in use in the clinic.

LITERATURE REFERENCES:

[1] Biology of incretins: GLP-1 and GIP; L.L. Baggio & D.J. Drucker; *Gastroenterology* **132**, 2131 (2007) • [2] GLP-1 (9-36) amide, cleavage product of GLP-1 (7-36) amide, is a glucoregulatory peptide; D. Elahi, et al.; *Obesity* **16**, 1501 (2008) • [3] Characterisation of the processing by human neutral endopeptidase 24.11 of GLP-1(7-36) amide and comparison of the substrate specificity of the enzyme for other glucagon-like peptides; K. Hupe-Sodmann, et al.; *Regul. Pept.* **58**, 149 (1995) • [4] Structural studies of a bifunctional inhibitor of neprilysin and DPP-IV; C. Oefner, et al.; *Acta Crystallogr. D Biol. Crystallogr.* **D63**, 975 (2007)

Glucagon

Glucagon

BML-GP1432-0100 100 µg

Human prepro-glucagon [51-81]. Blocking and immunizing peptide for the glucagon polyclonal antibody (BML-GA1181). **APPLICATIONS:** Standard for radioimmunoassay, coating peptide in enzyme immunoassays, complementary peptide for immunohistochemistry.

Glucagon, pAb

BML-GA1181-0025 25 µl

BML-GA1181-0100 100 µl

From rabbit. **IMMUNOGEN:** Natural pig pancreatic glucagons. **SPECIFICITY:** Recognizes mammalian glucagon. **APPLICATION:** IHC

Glucagon like Peptides [GLPs]

GLP-1, pAb

BML-GA1176-0025 25 µl

BML-GA1176-0100 100 µl

From rabbit. **IMMUNOGEN:** Synthetic peptide corresponding to aa 1-19 of human GLP-1. **SPECIFICITY:** Recognizes mammalian GLP1. **APPLICATION:** IHC. **BP:** BML-GP1436

GLP-2, pAb

BML-GA1179-0025 25 µl

BML-GA1179-0100 100 µl

From rabbit. **IMMUNOGEN:** Synthetic peptide corresponding to aa 1-33 of human GLP-2. **SPECIFICITY:** Recognizes mammalian GLP-2. **APPLICATION:** IHC. **BP:** BML-GP1439

GLP-1 (7-36) amide, pAb

BML-GA1178-0250 1 Vial

From rabbit. **IMMUNOGEN:** Synthetic peptide corresponding to aa 7-36 of human GLP-1. **SPECIFICITY:** Recognizes human, rat and other mammalian species GLP-1. **APPLICATION:** RIA. **BP:** BML-GP1438

DPPIV Proteins

DPPIV (human), (rec.)

BML-SE434-0025

25 mU

Active recombinant human DPPIV (aa 29-766). This represents a naturally-occurring cleaved (soluble) form of DPPIV. **BIOLOGICAL ACTIVITY:** Study of enzyme kinetics, cleave target substrates, and screen for inhibitors.

DPPIV (human), (rec.)

ALX-201-128-1

1 Vial

Produced in Sf9 cells.

DPPIV Kits

DPPIV assay kit for biological samples

BML-AK498-0001

1 Kit

The DPPIV assay kit for biological samples is a complete assay system designed to measure DPPIV activity in biological fluids such as plasma, serum, urine, and saliva. Uses for the kit include correlation of DPPIV activity with disease states or determination of the efficacy of DPPIV inhibitors administered *in vivo*. **QUANTITY:** 96 assays

DPPIV drug discovery kit

BML-AK499-0001

1 Kit

The DPPIV drug discovery kit is a complete assay system designed to screen inhibitors of DPPIV (DPP4; CD26). Using a convenient microplate format, DPPIV activity is monitored either colorimetrically, or fluorimetrically using substrates provided in the kit.

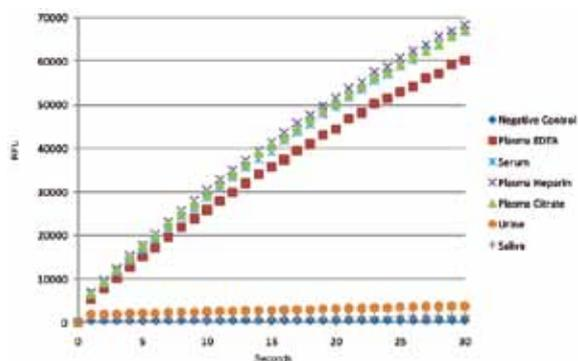


FIGURE: DPPIV activity in biological samples

HIGHLIGHT

DPP Substrates

GP-pNA

[H-Gly-Pro-p-nitroanilide]

BML-P188-0025

25 mg

Chromogenic substrate for dipeptidyl peptidase IV (DPIV, DPPIV, CD26), DPP7/II, DPP8, DPP9, FAP, PCP/PRCP, POP/PREP, and likely other prolyl peptidases. Can be used *in vitro* for inhibitor screening, on cultured cells, and in biological fluids. Hydrolysis of this substrate can be monitored at 405-410nm.

GP-AMC

[H-Gly-Pro-7-amino-4-methylcoumarin]

BML-P189-0005

5 mg

Fluorogenic substrate for dipeptidyl peptidase IV (DPIV, DPPIV, CD26) ($k_{cat}/K_m=2.6 \times 10^5 s^{-1} M^{-1}$), DPPII/VII, DPP8 and DPP9, FAP (fibroblast activation protein), and likely POP/PREP, but not glutamate carboxypeptidase II. Can be used *in vitro*, for inhibitor screening, and in biological fluids. Ex = 380nm, Em=460nm.

VP-AMC

BML-P448-0005

5 mg

Fluorogenic substrate likely cleaved by dipeptidyl peptidase IV (DPIV, DPPIV, CD26), DPP8, FAP (fibroblast activation protein), and other prolyl peptidases. Can be used *in vitro*, for inhibitor screening, and in biological fluids. Ex = 380 nm, Em = 460 nm.

DP-AMC

BML-P449-0005

5 mg

Fluorogenic substrate cleaved by dipeptidyl peptidase IV (DPIV, DPPIV, CD26) ($K_m=28\mu M$, $k_{cat}/K_m=3.7 \times 10^3 M^{-1} s^{-1}$), DPPII/VII ($K_m=1.12mM$, $k_{cat}/K_m=40.0 M^{-1} s^{-1}$), and likely DPP8 and other prolyl peptidases. Can be used *in vitro*, for inhibitor screening, and in biological fluids. Ex = 380 nm, Em = 460 nm.

WP-AMC

BML-P450-0005

5 mg

Fluorogenic substrate likely cleaved by dipeptidyl peptidase IV (DPIV, DPPIV, CD26), DPP8, and other prolyl peptidases. Can be used *in vitro*, for inhibitor screening, and in biological fluids. Ex = 380 nm, Em = 460 nm.

DPPIV [CD26] Antibody

DPPIV (human), pAb

BML-SA451-0100 100 µg

From rabbit. **IMMUNOGEN:** Synthetic peptide corresponding to a portion of human DPPIV. **SPECIFICITY:** Recognizes human DPPIV. **APPLICATION:** WB.

DPPIV Inhibitors

Diprotin A

[H-Ile-Pro-Ile-OH]

ALX-260-036-M005 5 mg

ALX-260-036-M025 25 mg

Synthetic. Inhibitor of dipeptidyl aminopeptidase IV.

LT: Diprotins A and B, inhibitors of dipeptidyl aminopeptidase IV, produced by bacteria: H. Umezawa, et al.; J. Antibiot. **37**, 422 (1984) • Purification and some properties of a membrane-bound dipeptidyl peptidase IV of guinea pig casein-induced intraperitoneal leukocytes: M. Kudo, et al.; J. Biochem. **97**, 1211 (1985) • Murine thymocytes possess specific cell surface-associated exoaminopeptidase activities: preferential expression by immature CD4- CD8- subpopulation: B. Bauvois; Eur. J. Immunol. **20**, 459 (1990)

Diprotin B

[H-Val-Pro-Leu-OH]

ALX-260-013-M005 5 mg

ALX-260-013-M025 25 mg

Synthetic. Inhibitor of dipeptidyl aminopeptidase IV.

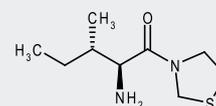
LT: Diprotins A and B, inhibitors of dipeptidyl aminopeptidase IV, produced by bacteria: H. Umezawa, et al.; J. Antibiot. **37**, 422 (1984)

P32/98

BML-PI142-0010 10 mg

BML-PI142-0050 50 mg

P32/98 (Ile-thiazolidide) is a specific, competitive transition-state substrate analog inhibitor of dipeptidyl peptidase IV (DPIV, DPPIV, CD26), with a K_i of 130nM. Useful in diabetes research. Has been used *in vivo* and in tissue culture.



P32/48

Related Products

FAP (human), (rec.)

BML-SE409-0010 10 µg

Active recombinant human FAP (aa 29-760), representing a naturally-occurring cleaved (soluble) form. **BIOLOGICAL ACTIVITY:** Study of enzyme kinetics, cleave target substrates, and screen for inhibitors.

POP/PREP (human), (rec.)

BML-SE545-0010 10 µg

Active recombinant human POP/PREP/PEP with an N-terminal purification tag. **BIOLOGICAL ACTIVITY:** Study of enzyme kinetics, cleave target substrates, and screen for inhibitors.

Nepriylsin

Nepriylsin (human), (rec.) (His-tag)

BML-SE532-0010 10 µg

Produced in Sf9 cells.

Nepriylsin (rat kidney), (purified)

BML-SE512-0010 10 µg

Isolated from rat kidney.

Mca-RPPGFSAFK(Dnp)

BML-P227-0001 1 mg

Fluorogenic substrate for ECE-1 and ECE-2, ACE1 and ACE2, neprilysin, MMP-2 and MMP-9, thimet oligopeptidase, as well as cathepsin A, cathepsin X/Z, BACE, and insulinase, but not MMP-1. Mca fluorescence is quenched by the Dnp group until cleavage separates them. Ex.: 328nm, Em.: 393nm, although the following Ex/Em have also been used: 320-340/400-420. This substrate is useful for inhibitor screening and kinetic analysis.

LT: Phosphinic tripeptides as dual angiotensin-converting enzyme C-domain and endothelin-converting enzyme-1 inhibitors: N. Jullien, et al.; J. Med. Chem. **53**, 208 (2010)

Nepriylsin (human), cAb (E19-P)

ALX-810-204-R100 100 µl

ALX-810-204-R500 500 µl

ALX-810-204-L001 1 ml

CLONE: E19-P. **ISOTYPE:** Rabbit IgG. **IMMUNOGEN:** Synthetic peptide corresponding to a C-terminal sequence of human neprilysin (CD10; CALLA). **SPECIFICITY:** Recognizes human neprilysin. **APPLICATION:** IHC (FS, PS).

Nepriylsin (human), cAb (E19-P) (Ready-to-Use)

ALX-810-404-L007 7 ml

ALX-810-404-L015 15 ml

CLONE: E19-P. **ISOTYPE:** Rabbit IgG. **IMMUNOGEN:** Synthetic peptide corresponding to a C-terminal sequence of human neprilysin. **SPECIFICITY:** Recognizes human neprilysin. **APPLICATION:** IHC (PS).

Antidiabetic Agents

Antidiabetic Agents

Genipin

ALX-350-383-M025 25 mg
ALX-350-383-M100 100 mg

Isolated from *Gardenia jasminoides Ellis*. Cell permeable inhibitor of uncoupling protein 2 (UCP2). Increases glucose-stimulated insulin secretion, mitochondrial membrane potential and ATP levels in pancreatic islet cells. Induces apoptosis. Protein cross-linking agent. Anti-inflammatory and anti-angiogenic.

LIT: Genipin-induced apoptosis in hepatoma cells is mediated by reactive oxygen species/c-Jun NH2-terminal kinase-dependent activation of mitochondrial pathway: B.C. Kim, et al.; *Biochem. Pharmacol.* **70**, 1398 (2005) • Genipin inhibits UCP2-mediated proton leak and acutely reverses obesity- and high glucose-induced beta cell dysfunction in isolated pancreatic islets: C.Y. Zhang, et al.; *Cell Metab.* **3**, 417 (2006) • Glucose sensing by POMC neurons regulates glucose homeostasis and is impaired in obesity: L.E. Parton, et al.; *Nature* **449**, 228 (2007)

Glyburide

[Glibenclamide]

BML-KC120-0001 1 g
BML-KC120-0005 5 g

An antidiabetic sulfonylurea which selectively blocks ATP-sensitive K⁺ channels. It blocks vascular smooth muscle relaxation produced by K_{ATP} channel openers. High affinity binding sites are widely distributed and have been identified in the cardiovascular system, the CNS and in pancreatic β-cells. It displays a large potency difference in pancreatic tissue (IC₅₀ in the nM range) versus vascular tissue (IC₅₀ in the μM range).

Glipizide

ALX-550-271-M100 100 mg
ALX-550-271-G001 1 g
ALX-550-271-G005 5 g

Antidiabetic agent. ATP-dependent K⁺ channel blocker. Blocks K⁺ channels in pancreatic β cells leading to an increase in calcium and insulin.

LIT: RP 49356 and cromakalim relax airway smooth muscle in vitro by opening a sulphonylurea-sensitive K⁺ channel: a comparison with nifedipine: D. Raeburn & T.J. Brown; *J. Pharmacol. Exp. Ther.* **256**, 480 (1991) • Quantitative autoradiography of the binding sites for [¹²⁵I] iodoglyburide, a novel high-affinity ligand for ATP-sensitive potassium channels in rat brain: D.R. Gehlert, et al.; *J. Pharmacol. Exp. Ther.* **257**, 901 (1991)

Metformin

ALX-270-432-G001 1 g
ALX-270-432-G005 5 g

Antidiabetic agent. Reduces the cardiovascular complications of diabetes. Does not induce hypoglycemia.

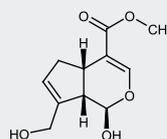
LIT: Effects of metformin and thiazolidinediones on suppression of hepatic glucose production and stimulation of glucose uptake in type 2 diabetes: a systematic review: A. Natali and E. Ferrannini; *Diabetologia* **49**, 434 (2006) • **For a comprehensive bibliography please visit our website.**

Pterostilbene

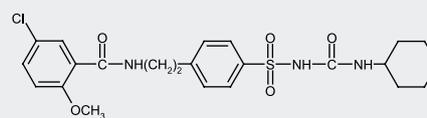
ALX-385-034-M025 25 mg
ALX-385-034-M100 100 mg
ALX-385-034-M500 500 mg

Synthetic. Cell permeable natural methoxylated analog of resveratrol (Prod. No. ALX-270-125). Antioxidant, antiproliferative, anti-inflammatory, anti-hyperglycemic and anti-diabetic agent. Induces apoptosis. Inhibits activator protein 1 (AP-1) and NF-κB activation. Moderately inhibits COX-1 and COX-2 (IC₅₀=19.8μM and 83.9μM).

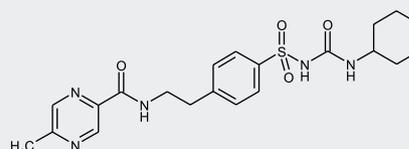
LIT: Pterostilbene, a new agonist for the peroxisome proliferator-activated receptor alpha-isoform, lowers plasma lipoproteins and cholesterol in hypercholesterolemic hamsters: A.M. Rimando, et al.; *J. Agric. Food Chem.* **53**, 3403 (2005) • Effect of pterostilbene on hepatic key enzymes of glucose metabolism in streptozotocin- and nicotinamide-induced diabetic rats: L. Pari & M.A. Satheesh; *Life Sci.* **79**, 641 (2006) • **For a comprehensive bibliography please visit our website.**



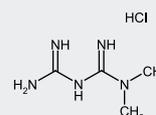
Genipin



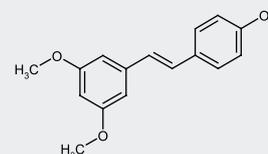
Glyburide



Glipizide



Metformin



Pterostilbene

HDACs

There is clear evidence of HDAC involvement in diabetes pathogenesis, and its pre-disposing factors, obesity and inflammation [1]. Development of adult type 2 diabetes in a rat model of intrauterine growth retardation is associated with HDAC1-induced epigenetic changes and these could be reversed by HDAC inhibition [2]. Due to its repression of GLUT4 expression in skeletal muscle, HDAC5 is implicated in regulation of blood glucose [3-5]. HDAC inhibitors have been found to prevent cytokine-induced toxicity in pancreatic β -cells [6] and to increase the pool of β -cells during development [7].

LITERATURE REFERENCES:

[1] Role of histone and transcription factor acetylation in diabetes pathogenesis: S.G. Gray & P. De Meys; *Diabetes Metab. Res. Rev.* **21**, 416 (2005) • [2] Development of type 2 diabetes following intrauterine growth retardation in rats is associated with progressive epigenetic silencing of Pdx1: J.H. Park, et al.; *J. Clin. Invest.* **118**, 2316 (2008) • [3] Restoration of insulin-sensitive glucose transporter (GLUT4) gene expression in muscle cells by the transcriptional coactivator PGC-1: L.F. Michael, et al.; *PNAS* **98**, 3820 (2001) • [4] Regulation of peroxisome proliferator-activated receptor gamma coactivator 1 alpha (PGC-1 alpha) and mitochondrial function by MEF2 and HDAC5: M.P. Czubyrt, et al.; *PNAS* **100**, 1711 (2003) • [5] AMP-activated protein kinase regulates GLUT4 transcription by phosphorylating histone deacetylase 5: S.L. McGee, et al.; *Diabetes* **57**, 860 (2008) • [6] Inhibition of histone deacetylases prevents cytokine-induced toxicity in beta cells: L. Larsen, et al. *Diabetologia* **50**, 779 (2007) • [7] Histone deacetylase inhibitors modify pancreatic cell fate determination and amplify endocrine progenitors: C. Haumaitre, et al.; *Mol. Cell Biol.* **28**, 6373 (2008)

HDAC Antibodies

Product	Specificity	Application	Prod. No.	Size
HDAC1, mAb (10E2)	Human and mouse	IHC, ICC, IP, WB	ALX-804-599-C200	200 μ g
HDAC1, pAb	Human, rat and mouse	IHC, IP, WB	BML-SA401-0100	100 μ g
HDAC2, pAb	Human, rat and mouse	IHC, IP, WB	BML-SA402-0100	100 μ g
HDAC3, pAb	Human, canine, hamster and rat	IP, WB	BML-SA403-0100	100 μ g
HDAC4, pAb	Human	WB	BML-SA404-0100	100 μ g
HDAC4 (NT), pAb	Human and mouse	WB	ALX-210-339-C100	100 μ g
HDAC5, pAb	Human and mouse	IP, WB	ALX-210-340-C100	100 μ g
HDAC6, pAb	Human and mouse	IP, WB	ALX-210-341-C100	100 μ g

HDAC Inhibitors

Trichostatin A

BML-GR309-0001 1 mg
BML-GR309-0005 5 mg

Trichostatin A (TSA) is a potent and reversible inhibitor of histone deacetylases. In HeLa cells, TSA blocked cell cycle progression at G1 and induced a 12-fold increase in intracellular levels of gelsolin. In cells latently infected with HIV-1, TSA induced the transcriptional activation of the HIV-1 promoter, which resulted in a marked increase in virus production. In NIH 3T3 cells, TSA induced reversion of oncogenic ras-transformed cells to a normal morphology. In Jurkat cells, TSA inhibited IL-2 gene expression (IC_{50} = 73nM) and displayed immunosuppressive activity in a mouse model. Induces increased acetylation of GATA4, a cardiac-specific transcription factor and increases cardiac muscle cell differentiation. Trichostatin A is a useful tool for induction of hyperacetylation of cellular histones and for further elucidation of their role in gene expression.

LIT: Histone deacetylase: a regulator of transcription: A.P. Wolffe; *Science* **272**, 371 (1996) • Trichostatin A induces morphological changes and gelsolin expression by inhibiting histone deacetylase in human carcinoma cell lines: Y. Hoshikawa, et al.; *Exp. Cell Res.* **214**, 189 (1994) • Trichostatin A inhibits both ras-induced neurite outgrowth of PC12 cells and morphological transformation of NIH3T3 cells: M. Futamura, et al.; *Oncogene* **10**, 1119 (1995) • Selective inhibition of IL-2 gene expression by trichostatin A, a potent inhibitor of mammalian histone deacetylase: I. Takahashi, et al.; *J. Antibiot. (Tokyo)* **49**, 453 (1996)

Suberoyl bis-hydroxamic acid

BML-GR323-0100 100 mg
BML-GR323-0500 500 mg

Induces cellular differentiation. Inhibits histone deacetylases.

LIT: Cytodifferentiating agents affect the replication of herpes simplex virus type 1 in the absence of functional VP16: C.M. Preston & McFarlane, M.; *Virology* **249**, 418 (1998)

Depudecin

BML-EI319-0100 100 μ g

Depudecin is a fungal metabolite that reverts ras- and src-transformed NIH3T3 cells to a flat phenotype (1 μ g/ml). It is a potent HDAC inhibitor (IC_{50} = 4.7 μ M for HDAC1) and displays anti-angiogenic activity.

LIT: Depudecin, a microbial metabolite containing two epoxide groups, exhibits anti-angiogenic activity in vivo: T. Oikawa et al.; *Biol. Pharm. Bull.* **18**, 1305 (1995) • Depudecin induces morphological reversion of transformed fibroblasts via the inhibition of histone deacetylase: H.J. Kwon et al.; *PNAS* **95**, 3356 (1998) • Synthesis and cellular characterization of the detransformation agent, (-)-depudecin: J. Shimada et al.; *Chem. Biol.* **2**, 517 (1995) • Depudecin: a novel compound inducing the flat phenotype of NIH3T3 cells doubly transformed by ras- and src-oncogene, produced by *Alternaria brassicicola*: M. Matsumoto et al.; *J. Antibiot. (Tokyo)* **45**, 879 (1992)

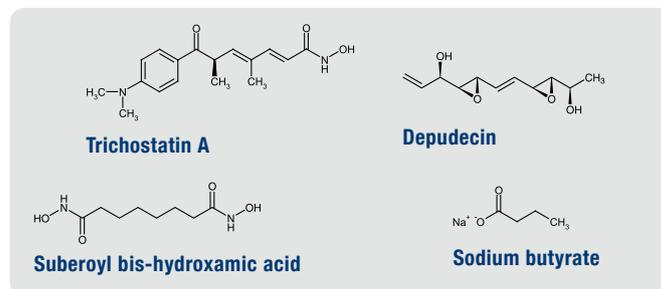
Sodium butyrate

[Butyric acid . sodium salt]

ALX-270-301-G001 1 g

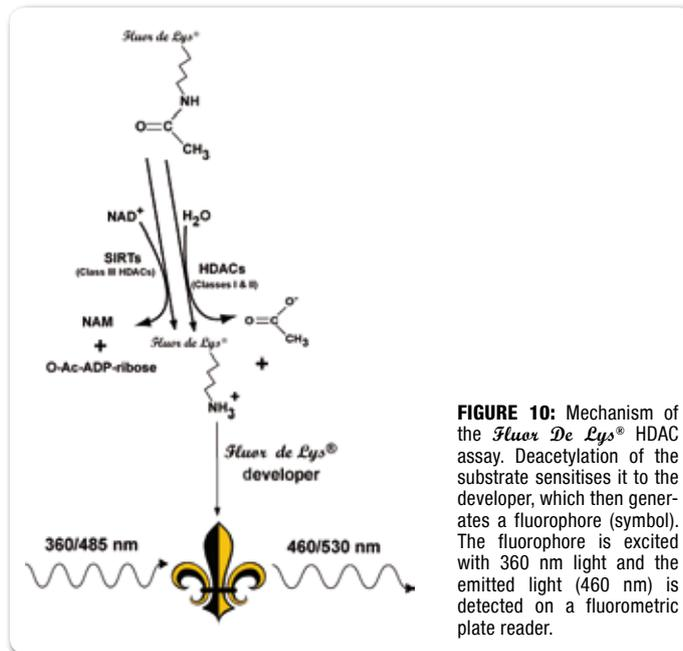
Short-chain fatty acid shown to be an inhibitor of histone deacetylases (HDACs). Induces apoptosis.

LIT: Sodium butyrate inhibits histone deacetylation in cultured cells: E.P. Candido, et al.; *Cell* **14**, 105 (1978) • Reduction of telomerase activity in human liver cancer cells by a histone deacetylase inhibitor: M. Nakamura, et al.; *J. Cell Physiol.* **187**, 392 (2001) • Effects of retinoic acid and sodium butyrate on gene expression, histone acetylation and inhibition of proliferation of melanoma cells: K. Demary, et al.; *Cancer Lett.* **163**, 103 (2001) • **For a comprehensive bibliography please visit our website.**



Activity Assay Kits

The *Fluor de Lys*[®] (Fluorescent deacetylation of lysine) Assay System was developed for simple, nonradioactive measurement of deacetylase activity. Deacetylation of a *Fluor de Lys*[®] substrate by a source of deacetylase activity, including purified enzyme or enzyme complex, cell lysate or whole cells, sensitizes the substrate so that, in a second step, treatment with the appropriate *Fluor de Lys*[®] Developer produces a fluorophore (see Figure 10). The assay is compatible with class I, IIb & IV HDACs and sirtuins and a variety of substrates are available based on acetylated sites found in p53 and histones. The new *Fluor de Lys*[®]-Green substrate (incorporated into BML-AK530) offers a higher sensitivity and an excitation and emission (485/530nm) that avoids quenching and fluorescent interference from compounds absorbing in the near UV and blue range. *Fluor de Lys*[®] Substrate is cell-permeable and allows cell-based determination of HDAC activity (Prod. No. BML-AK503). The *Fluor de Lys*[®] system is the basis of several activity assays/drug discovery kits for HDACs (Prod. No.s BML-AK500, BML-AK511, BML-AK516, BML-AK518, BML-AK530 & BML-AK531) and sirtuins (BML-AK555, BML-AK556 and BML-AK557) that allows straightforward screening for new modulators of deacetylase activity.



Selected citations using the *Fluor de Lys*[®] activity assays:

X. Zhou et al.; PNAS **98**, 10572 (2001) • K.J. Bitterman et al.; J. Biol. Chem. **277**, 45099 (2002) • B. Heltweg and Jung, M.; Anal. Biochem. **302**, 175 (2002) • K. Ito et al.; PNAS **99**, 8921 (2002) • S. Milutinovic et al.; J. Biol. Chem. **277**, 20974 (2002) • R.M. Anderson et al.; Science **302**, 2124 (2003) • K.T. Howitz et al.; Nature **425**, 191 (2003) • G.V. Kapustin et al.; Org. Lett. **5**, 3053 (2003) • D.K. Kim et al.; J. Med. Chem. **46**, 5745 (2003) • C.G. Kleer et al.; PNAS **100**, 11606 (2003) • K. Zhao et al.; Nat. Struct. Biol. **10**, 864 (2003) • T. Suzuki et al.; Bioorg. Med. Chem. Lett. **13**, 4321 (2003) • B.G. Cosio et al.; Am. J. Respir. Crit Care Med. **170**, 141 (2004) • C.M. Gallo et al.; Mol. Cell Biol. **24**, 1301 (2004) • N. Gurvich et al.; Cancer Res. **64**, 1079 (2004) • L.H. Wang et al.; Nat. Med. **10**, 40 (2004) • J.G. Wood et al.; Nature **430**, 686 (2004) • F. Yeung et al.; EMBO J. **23**, 2369 (2004) • J.L. Avalos et al.; Mol. Cell **17**, 855 (2005) • K. Ito et al.; N. Engl. J. Med. **352**, 1967 (2005) • A. Mai et al.; J. Med. Chem. **48**, 7789 (2005) • E. Michishita et al.; Mol. Biol. Cell **16**, 4623 (2005) • A.D. Napper et al.; J. Med. Chem. **48**, 8045 (2005) • T. Suzuki et al.; J. Med. Chem. **48**, 1019 (2005) • P. Aksoy et al.; BBRC **349**, 353 (2006) • V.C. de Boer et al.; Mech. Ageing Dev. **127**, 618 (2006) • S.L. Gantt et al.; Biochemistry **45**, 6170 (2006) • W. Gu et al.; Bioorg. Med. Chem. **14**, 3320 (2006) • D. Herman et al.; Nat. Chem. Biol. **2**, 551 (2006) • X. Li et al.; Cancer Res. **66**, 9323 (2006) • V.M. Nayagam et al.; J. Biomol. Screen. **11**, 959 (2006) • J.M. Solomon et al.; Mol. Cell Biol. **26**, 28 (2006) • P.H. Kiviranta et al.; Bioorg. Med. Chem. Lett. **17**, 2448 (2007) • T.F. Outeiro et al.; Science **317**, 516 (2007) • S. Lain et al.; Cancer Cell **13**, 454 (2008) • B. Jung-Hynes et al.; J. Biol. Chem. **284**, 3823 (2009)

Product	Prod. No.	Size
<i>Fluor-de-Lys</i> [®] HDAC fluorometric activity assay kit	BML-AK500-0001	1 Kit
Color-de-Lys [™] HDAC colorimetric activity assay kit	BML-AK501-0001	1 Kit
<i>Fluor-de-Lys</i> [®] HDAC1 fluorometric drug discovery assay kit	BML-AK511-0001	1 Kit
<i>Fluor-de-Lys</i> [®] HDAC8 fluorometric drug discovery kit	BML-AK518-0001	1 Kit
<i>Fluor-de-Lys</i> [®] HDAC fluorometric cellular activity assay kit	BML-AK503-0001	1 Kit
<i>Fluor-de-Lys</i> [®] HDAC3/NCOR1 fluorometric drug discovery kit	BML-AK531-0001	1 Kit
<i>Fluor-de-Lys</i> [®] -Green HDAC2 fluorometric drug discovery assay kit	BML-AK512-0001	1 Kit
<i>Fluor-de-Lys</i> [®] -Green HDAC fluorometric activity assay kit	BML-AK530-0001	1 Kit
<i>Fluor-de-Lys</i> [®] HDAC6 fluorometric drug discovery kit	BML-AK516-0001	1 Kit

Sirtuins [SIRT1s]

Sirtuins perform multiple regulatory functions in gene expression and metabolism via their action on histones, transcription factors, structural proteins and enzymes [1]. An enzyme family consisting mostly of NAD⁺-dependent lysine deacetylases (class III HDACs) [2], the sirtuins also include some ADP-ribosyl transferases [3]. All sirtuins use one important metabolic cofactor, NAD⁺, as a cosubstrate and the acetylation/deacetylation cycle occurs at the expense of another, acetyl-CoA. It has therefore been suggested that sirtuins sense the cell's energetic state, for example via NAD⁺ and/or nicotinamide concentrations [4, 5] or the NAD⁺:NADH ratio [6], and that their activity is regulated accordingly. Several of the seven members of the human sirtuin family (SIRT1-SIRT7) [2] are known to be profoundly involved in the regulation of insulin signaling [3, 7-9] and energy metabolism [10-18] and are therefore considered promising therapeutic targets for diabetes and obesity [19]. In rodent studies, SIRT1 activators such as resveratrol [20] and SIRT1720 [21] have been shown to improve insulin sensitivity [17, 21-23], lower plasma glucose [17, 21-23], increase mitochondrial function [17, 21-23] and muscle aerobic capacity [17, 23] and to protect against diet-induced obesity and/or its associated pathologies [17, 22, 23].

LITERATURE REFERENCES:

[1] Sirtuins in mammals: insights into their biological function: S. Michan & D. Sinclair; *Biochem. J.* **404**, 1 (2007) • [2] Phylogenetic classification of prokaryotic and eukaryotic Sir2-like proteins: R.A. Frye; *BBRC* **273**, 793 (2000) • [3] SIRT4 inhibits glutamate dehydrogenase and opposes the effects of calorie restriction in pancreatic beta cells: M.C. Haigis, et al.; *Cell* **126**, 941 (2006) • [4] Requirement of NAD and SIRT2 for life-span extension by calorie restriction in *Saccharomyces cerevisiae*: S.J. Lin, et al.; *Science* **289**, 2126 (2000) • [5] Yeast life-span extension by calorie restriction is independent of NAD fluctuation: R.M. Anderson, et al.; *Science* **302**, 2124 (2003) • [6] Calorie restriction extends yeast life span by lowering the level of NADH: S.J. Lin, et al.; *Genes Dev.* **18**, 12 (2004) • [7] Sirt1 regulates insulin secretion by repressing UCP2 in pancreatic beta cells: L. Bordone, et al.; *PLoS Biol.* **4**, e31 (2006) • [8] SIRT1 improves insulin sensitivity under insulin-resistant conditions by repressing PTP1B: C. Sun, et al.; *Cell Metab.* **6**, 307 (2007) • [9] The direct involvement of SirT1 in insulin-induced insulin receptor substrate-2 tyrosine phosphorylation: J. Zhang; *J. Biol. Chem.* **282**, 34356 (2007) • [10] Sirt1 promotes fat mobilization in white adipocytes by repressing PPAR-gamma: F. Picard, et al.; *Nature* **429**, 771 (2004) • [11] Sirtuins deacetylate and activate mammalian acetyl-CoA synthetases: W.C. Hallows, et al.; *PNAS* **103**, 10230 (2006) • [12] Reversible lysine acetylation controls the activity of the mitochondrial enzyme acetyl-CoA synthetase 2: B. Schwer, et al.; *PNAS* **103**, 10224 (2006) • [13] SIRT2 regulates adipocyte differentiation through FoxO1 acetylation/deacetylation: E. Jing, et al.; *Cell Metab.* **6**, 105 (2007) • [14] Calorie restriction promotes mammalian cell survival by inducing the SIRT1 deacetylase: H.Y. Cohen, et al.; *Science* **305**, 390 (2004) • [15] Tissue-specific regulation of SIRT1 by calorie restriction: D. Chen, et al.; *Genes Dev.* **22**, 1753 (2008) • [16] Conserved metabolic regulatory functions of sirtuins: B. Schwer & E. Verdin; *Cell Metab.* **7**, 104 (2008) • [17] Resveratrol improves mitochondrial function and protects against metabolic disease by activating SIRT1 and PGC-1alpha: M. Lagouge, et al.; *Cell* **127**, 1109 (2006) • [18] Regulation of SIRT1 protein levels by nutrient availability: Y. Kanfi, et al.; *FEBS Lett.* **582**, 2417 (2008) • [19] Sirtuins--novel therapeutic targets to treat age-associated diseases: S. Lavu, et al.; *Nat. Rev. Drug Discov.* **7**, 841 (2008) • [20] Small molecule activators of sirtuins extend *Saccharomyces cerevisiae* lifespan: K.T. Howitz, et al.; *Nature* **425**, 191 (2003) • [21] Small molecule activators of SIRT1 as therapeutics for the treatment of type 2 diabetes: J.C. Milne, et al.; *Nature* **450**, 712 (2007) • [22] Resveratrol improves health and survival of mice on a high-calorie diet: J.A. Baur, et al.; *Nature* **444**, 337 (2006) • [23] Specific SIRT1 activation mimics low energy levels and protects against diet-induced metabolic disorders by enhancing fat oxidation: J.N. Feige, et al.; *Cell Metab.* **8**, 347 (2008)

Sirtuin Kits

Fluor-de-Lys[®] SIRT1 fluorometric drug discovery assay kit

BML-AK555-0001

1 Kit

The SIRT1 fluorescent activity assay is based on the unique *Fluor de Lys*[®]-SIRT1 substrate/developer II combination. The *Fluor de Lys*[®]-SIRT1 substrate is a peptide comprising amino acids 379-382 of human p53 (Arg-His-Lys-Lys(Ac)). The assay's fluorescence signal is generated in proportion to the amount of deacetylation of the lysine corresponding to Lys³⁸², a known *in vivo* target of SIRT1 activity. *Fluor de Lys*[®]-SIRT1 was the substrate deacetylated most efficiently by SIRT1 from among a panel of substrates patterned on p53, histone H3 and histone H4 acetylation.

Fluor-de-Lys[®] SIRT2 fluorometric drug discovery assay kit

BML-AK556-0001

1 Kit

The SIRT2 fluorescent activity assay is based on the unique *Fluor de Lys*[®] substrate/developer II combination. The *Fluor de Lys*[®]-SIRT2 substrate is unique peptide based on amino acids 317-320 of p53 (Gln-Pro-Lys-Lys(Ac)), a site of regulatory acetylation by PCAF acetyltransferase (Lys³²⁰). It was selected as the best substrate for SIRT2 from a panel of substrates patterned on p53, histone H3, and histone H4.

Fluor-de-Lys[®] SIRT3 fluorometric drug discovery assay kit

BML-AK557-0001

1 Kit

The SIRT3 fluorimetric assay system is based on the *Fluor de Lys*[®] substrate/developer II combination. The *Fluor de Lys*[®]-SIRT2 substrate is unique peptide based on amino acids 317-320 of p53 (Gln-Pro-Lys-Lys(Ac)), a site of regulatory acetylation by PCAF acetyltransferase (Lys³²⁰). It was selected as the best substrate for SIRT3 from a panel of substrates patterned on p53, histone H3, and histone H4.

Visit www.enzolifesciences.com for a complete offering of sirtuin enzymes, substrates and inhibitors.

Sirtuin Antibodies

Product	Specificity	Application	Prod. No.	Size
SIRT1 (human), pAb	Human	WB	BML-SA427-0100	100 µl
SIRT2 (human), pAb	Human	WB	BML-SA444-0100	100 µl
SIRT3, pAb	Human, rat and bovine	WB	BML-SA463-0100	100 µl
SIRT5, pAb	Human, mouse, rat and bovine	WB	BML-SA464-0100	100 µl

Sirtuin Activators

Butein

ALX-350-246-M010 10 mg

Plant polyphenol. Activator of human deacetylase SIRT1. Specific tyrosine kinase inhibitor. Potently inhibits the tyrosine kinase activity of the EGF receptor and p60c-src. Potent antioxidant and anti-inflammatory agent. Inhibits glutathione reductase and rat liver glutathione S-transferase. Inhibits aromatase, showing chemopreventive properties. Directly inhibits IKK.

Lit: Butein, a specific protein tyrosine kinase inhibitor: E.-B. Yang, et al.; *BBRC* **245**, 435 (1998) • Flavonoids inhibit cell growth and induce apoptosis in B16 melanoma 4A5 cells: K. Iwashita, et al.; *Biosci. Biotechnol. Biochem.* **64**, 1813 (2000) • Small molecule activators of sirtuins extend *Saccharomyces cerevisiae* lifespan: K.T. Howitz, et al.; *Nature* **425**, 191 (2003) • The plant polyphenol butein inhibits testosterone-induced proliferation in breast cancer cells expressing aromatase: Y. Wang, et al.; *Life Sci.* **77**, 39 (2005) • Butein, a tetrahydroxychalcone, inhibits nuclear factor (NF)-kappaB and NF-kappaB-regulated gene expression through direct inhibition of I kappaBalpha kinase beta on cysteine 179 residue: M.K. Pandey, et al.; *J. Biol. Chem.* **282**, 17340 (2007) • **For a comprehensive bibliography please visit our website.**

Piceatannol

ALX-270-202-M001 1 mg

ALX-270-202-M005 5 mg

ALX-270-202-M010 10 mg

ALX-270-202-M050 50 mg

Synthetic. Activator of human deacetylase SIRT1. Originally isolated from *Euphorbia lagascae*. Selective protein tyrosine kinase Syk inhibitor.

Lit: Piceatannol, a stilbene phytochemical, inhibits mitochondrial F₀F₁-ATPase activity by targeting the F₁ complex: J. Zheng & V.D. Ramirez; *BBRC* **261**, 499 (1999) • Microarray analysis of piceatannol-induced changes in gene expression in human gastric cancer cells: D. Jeoung, et al.; *Biotechnology Lett.* **24**, 463 (2002) • Small molecule activators of sirtuins extend *Saccharomyces cerevisiae* lifespan: K.T. Howitz, et al.; *Nature* **425**, 191 (2003) • **For a comprehensive bibliography please visit our website.**

Quercetin . dihydrate

ALX-385-001-G005 5 g

ALX-385-001-G025 25 g

Isolated from *Sophora japonica* L. Activator of human deacetylase SIRT1. Antioxidant flavonoid. Inhibitor of mitochondrial ATPase, cAMP- and cGMP-phosphodiesterases. Inhibitor of protein tyrosine kinases and protein kinase C (PKC). Induces apoptosis. Blocks cells at the G₀/G₁ interface. Reversible inhibitor of fatty acid synthase (FAS). Inhibits the production of the inflammatory mediators nitric oxide (NO), TNF- α and IL-12 in activated macrophages

Lit: Induction of apoptosis by quercetin: involvement of heat shock protein: Y.Q. Wei, et al.; *Cancer Res.* **54**, 4952 (1994) • Molecular mechanisms in the antiproliferative action of quercetin: B. Csokay, et al.; *Life Sci.* **60**, 2157 (1997) • Small molecule activators of sirtuins extend *Saccharomyces cerevisiae* lifespan: K.T. Howitz, et al.; *Nature* **425**, 191 (2003) • Pharmacological inhibitors of Fatty Acid Synthase (FASN)--catalyzed endogenous fatty acid biogenesis: a new family of anticancer agents?: R. Lupu & J. A. Menendez; *Curr. Pharm. Biotechnol.* **7**, 483 (2006) • **For a comprehensive bibliography please visit our website.**

Resveratrol

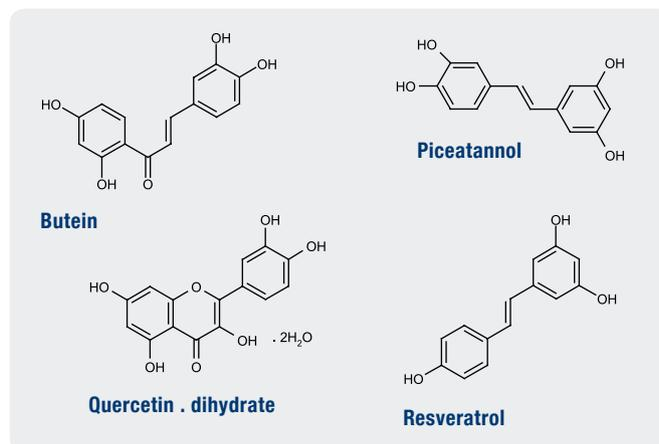
[*trans*-3,4',5-Trihydroxystilbene]

BML-FR104-0100 100 mg

BML-FR104-0500 500 mg

Phenolic antioxidant found in grapes and wine. Potent activator of human deacetylase SIRT1 (maximum effect at 100-200 μ M). Improves insulin sensitivity in rodent models. Increases longevity in yeast. Displays antiproliferative effect on human breast epithelial cells. Displays a variety of pharmacological activities including: estrogen receptor agonist activity, inhibition of ribonucleotide reductase and DNA synthesis, inhibition of eicosanoid synthesis, cell cycle arrest at S/G₂ phase and induction of apoptosis. Inhibits PKD in vitro (IC₅₀=200 μ M) and *in vivo* (IC₅₀=800 μ M), but not PKC isoforms. Protects against 4-hydroxynonenal (4-HNE) induced oxidative stress and apoptosis. Shows cancer chemopreventive activity. Specific inhibitor of cyclooxygenase-1 (COX-1). Inhibits the hydroperoxidase activity of COX-1.

Lit: Cancer chemopreventive activity of resveratrol, a natural product derived from grapes: M. Jang, et al.; *Science* **275**, 218 (1997) • Small molecule activators of sirtuins extend *Saccharomyces cerevisiae* lifespan: K.T. Howitz, et al.; *Nature* **425**, 191 (2003) • Resveratrol improves mitochondrial function and protects against metabolic disease by activating SIRT1 and PGC-1 α : M. Lagouge, et al.; *Cell* **127**, 1109 (2006) • Small molecule activators of SIRT1 as therapeutics for the treatment of type 2 diabetes: J.C. Milne, et al.; *Nature* **450**, 712 (2007) • Resveratrol improves health and survival of mice on a high-calorie diet: J.A. Baur, et al.; *Nature* **444**, 337 (2006) • Specific SIRT1 activation mimics low energy levels and protects against diet-induced metabolic disorders by enhancing fat oxidation: J.N. Feige, et al.; *Cell Metab.* **8**, 347 (2008) • Resveratrol interferes with AKT activity and triggers apoptosis in human uterine cancer cells: E. Sexton, et al.; *Mol. Cancer* **5**, 45 (2006) • Resveratrol and its analogs: Defense against cancer, coronary disease and neurodegenerative maladies or just a fad?: P. Saiko, et al.; *Mutat. Res.* **658**, 68 (2008) • **For a comprehensive bibliography please visit our website.**



Epigenetics & Chromatin Modification Catalog

Features a comprehensive range of over 150 products, including histone deacetylases (HDACs) and sirtuins (enzymes, substrates, activators, inhibitors, and antibodies).

Visit www.enzolifesciences.com for a complete listing or ask for a free copy of our new product flyers.



International Distributors

Argentina

LAB SCIENTIFIC, INC. (USA)
Tel: (305) 716-9922
Fax: (305) 716-9923
E-mail: labscientific@labscent.com

Australia

SAPPHIRE BIOSCIENCE Pty. Ltd.
Tel: +61 (0) 2 9698 2022
Fax: +61 (0) 2 9698 1022
E-mail: sales@sapphirebioscience.com

UNITED BIORESEARCH PRODUCTS
Tel: +61 (0) 2 9651 3736
Fax: +61 (0) 2 9651 4247
E-mail: kirrily@unitedbioresearch.com.au

Austria

EUBIO
Tel: (0)1 89 50 145
Fax: (0)1 89 50 145-14
E-mail: koeck@eubio.at

Bangladesh

FUTURE BUSINESS VISION
Tel: (0)2 8631173
Fax: (0)2 8651847
E-mail: salam.fbv@gmail.com

Belarus

CHIMMED Inc.
Tel: +7 095 728 4192
Fax: +7 095 742 8341
E-mail: bio@chimmed.ru

Belgium

ENZO LIFE SCIENCES BVBA
Tel: +32 (0) 3 466 04 20
Fax: +32 (0) 3 466 04 29
E-mail: info-be@enzolifesciences.com

Bosnia & Herzegovina

A-Ž LAB D.O.O.
Tel: +386 (0)1 433 63 22 /
+386 (0)1 230 18 84
Fax: +386 (0)1 230 19 85
E-mail: az.consulting@siol.net

Brazil

LGC BIOTECHNOLOGIA
Tel: (0)21 2583 4268 / (0)21 3273 5828
Fax: (0)21 3273 5896
E-mail: info@lgcbio.com.br

SELLEX S.A.C.
Tel: (0)11 5506 4646
Fax: (0)11 5507 4204
E-mail: vendas@sellex.com

Canada

ENZO LIFE SCIENCES INTERNATIONAL, INC.
Tel: (610) 941-0430
Toll Free Tel: 1-800-942-0430
Fax: (610) 941-9252
E-mail: info-usa@enzolifesciences.com

CEDARLANE LABORATORIES

Tel: (289) 288-0001
Toll Free: 1-800-268-5058
Fax: (289) 288-0020
Toll Free: 1-800-638-5099
E-mail: general@cedarlanelabs.com

MJS BioLynx Inc.

Tel: (613) 498-2126
Toll Free: 1-888-593-5969
Fax: (613) 342-1341
E-mail: tech@biolynx.ca

Chile

BIOCANT Ltda.
Tel: (0)2 683 2437
Fax: (0)56 2 683 8823
E-mail: info@biocant.cl

China

BEIJING BITAB BIOTECH Co. Ltd.
Tel: (0)10 8201 5225
Fax: (0)10 6201 5131
E-mail: info@bitebo.com

BOPPARD
Tel: (0)21 6288 4751 (SH)
Tel: (0)20 8732 6381 (GZ)
Fax: (0)21 6288 4752 (SH)
Fax: (0)20 8732 6382 (GZ)
E-mail: info@boppard.cn

GENETIMES TECHNOLOGY Inc.
Tel: (0)21 5426 2677
Fax: (0)21 6439 8855
E-mail: order@genetimes.com.cn

ITS CHINA
Tel: (0)21 648 144 28/98
Fax: (0)21 643 93402
E-mail: info@its-science-china.com

KANGCHEN BIO-TECH
Tel: (0)21 6455 1989
Toll Free: 800 820 5058 (China only)
Fax: (0)21 6455 2021
E-mail: order@kangchen.com.cn

MULTISCIENCES BIOTECH Co. Ltd.
Tel: (0)571 8816 3301
Toll Free: (0)800 8571 184
Fax: (0)571 8816 3303
E-mail: service@gotofcm.com

NEOBIO SCIENCE TECHNOLOGY
Tel: (0)755 26 755 677
Hot line: 4006 800 892
Fax: (0)755 26 755 877
E-mail: info@neobioscience.com

TWC BIOSEARCH INTERNATIONAL
Tel: (0)852 2649 9988
Fax: (0)852 2635 0379
E-mail: support@twcbiosearch.com

Colombia

LAB SCIENTIFIC, INC. (USA)
Tel: (305) 716-9922
Fax: (305) 716-9923
E-mail: labscientific@labscent.com

Cyprus

SB BIOTECHNOLOGY SUPPLIERS SA
Tel: +30 210 823 3373 / +30 210 691 0148
Fax: +30 210 825 9987
E-mail: info@sbbiotech.gr

Czech Republic

GENETICA s.r.o.
Tel: +420 2 7270 1055
Fax: +420 2 7270 1739
E-mail: geneticka@genetica.cz

Denmark

SMS GRUPPEN
Tel: (0)4586 4400
Fax: (0)4586 4881
E-mail: mail@sms-gruppen.dk

Ecuador

LAB SCIENTIFIC, INC. (USA)
Tel: (305) 716-9922
Fax: (305) 716-9923
E-mail: rshlesinger@labscent.com

Egypt

NEW TEST Co. (NTCo)
Tel: (0)3544 4736
Fax: (0)3359 6836
E-mail: info@newtest.com.eg

Estonia

IN VITRO EESTI OÜ
Tel: +372 630 65 20
Fax: +372 630 65 22
E-mail: info@invitro.ee

Finland

NUPPULINNA LABORATORIOPALVELU OY
Tel: (0)20 792 0350
Fax: (0)20 792 0351
E-mail: nuppulinna@dlc.fi

France

ENZO LIFE SCIENCES FRANCE
c/o Covalab s.a.s.
Tel: +33/0 472 440 655
Fax: +33/0 437 484 239
E-mail: info-fr@enzolifesciences.com

Germany

ENZO LIFE SCIENCES GmbH
Tel: (0)7621 5500 522
Toll Free 0800 253 9472
Fax: (0)7621 5500 527
E-mail: info-de@enzolifesciences.com

Greece

SB BIOTECHNOLOGY SUPPLIERS SA
Tel: +30 210 823 3373 / +30 210 691 0148
Fax: +30 210 825 9987
E-mail: info@sbbiotech.gr

Hong Kong

BOPPARD (H.K) Co. Ltd
Tel: +852 2799 9019
Fax: +852 2799 9808
E-mail: info@boppard.com.hk

Hungary

BIOMARKER Ltd
Tel: (0)28 419 986
Fax: (0)28 422 319
E-mail: biomarker@biomarker.hu

India

HYSEL INDIA Pvt. Ltd.
Tel: (0)11 2622 7801/02/03/04
Fax: (0)11 2622 7805
E-mail: hysel@del2.vsnl.net.in

GAURAV ENTERPRISE (AGRA)
Tel: (0)562 288 3724
Fax: (0)562 288 1414
E-mail: girish640@yahoo.co.in

LABEX CORPORATION
Tel: (0)11 2612 4727 / (0)11 2613 5922 /
(0)11 41771988
Fax: (0)11 2612 4735 / (0)11 2689 3172
E-mail: labex@labex.net

PRO LAB MARKETING Pvt. Ltd.
Tel: (0)11 6660 7725 / (0)11 6565 2166
Fax: (0)11 6660 7726 / (0)11 4165 8854
E-mail: info@prolabmarketing.com

Indonesia

ITS INDONESIA
Tel: (0)21 451 6222
Fax: (0)21 451 6223
E-mail: info@its-indonesia.com

Iran

HORMOZ PAJOHAN LAB. EQUIPMENT Ltd.
Tel: (0)21 8888 3444
Fax: (0)21 8877 0192
E-mail: ahmadi@hermes-pajohan.com

Iraq

IRAQ HEART Co. Ltd.
Tel: (0)790 171 7504
Cell: (0)770 278 7372 / (0)780 780 9800
E-mail: maithem_ghsan@yahoo.com

Ireland

ENZO LIFE SCIENCES (UK) LTD.
Tel: 0845 601 1488 / +44/0 1392 825900
Fax: +44/0 1392 825910
E-mail: info-uk@enzolifesciences.com

Israel

ALMOG DIAGNOSTIC & MEDICAL EQUIPMENT Ltd.
Tel: (0)3977 3390
Fax: (0)3977 3391
E-mail: info@almog.co.il

GADOT LABORATORY SUPPLY Ltd.
Tel: (0)5075 222 49
Toll Free: 1 800 20 22 20 (Israel Only)
Fax: 1 800 300 707
E-mail: mira@gadot.com

Italy

VINCI-BIOCHEM
Tel: (0)571 568 147
Fax: (0)571 568 132
E-mail: vb@vincibiochem.it

Japan

BIOLINKS K.K.
Tel: (0)3 5443 6891
Fax: (0)3 5443 0271
E-mail: info@biolinks.co.jp

COSMO BIO Co. Ltd.
Tel: (0)3 5632 9610
Fax: (0)3 5632 9619
E-mail: mail@cosmobio.co.jp

FUNAKOSHI Co., Ltd.
Tel: (0)3 5684 1620
Fax: (0)3 5684 1775
E-mail: reagent@funakoshi.co.jp

Kazakhstan

CHIMMED Inc.
Tel: +7 095 728 4192
Fax: +7 095 742 8341
E-mail: bio@chimmed.ru

Korea, South

CHUN YANG TECH
Tel: +82 32 624 0160 2
Fax: +82 32 624 0163
E-mail: 123ky@naver.com

SERVLAB CO.
Tel: +82 2 449 8787
Fax: +82 2 499 8786
E-mail: servlab@servlab.co.kr

Latvia

IN VITRO EESTI OÜ
Tel: +372 630 65 20
Fax: +372 630 65 22
E-mail: info@invitro.ee

Lithuania

IN VITRO EESTI OÜ
Tel: +372 630 65 20
Fax: +372 630 65 22
E-mail: info@invitro.ee

Luxembourg

ENZO LIFE SCIENCES BVBA
Tel: +32 (0) 3 466 04 20
Fax: +32 (0) 3 466 04 29
E-mail: info-be@enzolifesciences.com

Malaysia

INTERSCIENCE SDN BHD
Tel: (0)3 5740 9888
Fax: (0)3 5740 9866
E-mail: info@its-interscience.com

Mexico

CONSULTORIA DE LABORATORIOS S.A.
Tel: +52 (0) 55 4622 2691
Fax: +52 (0) 55 4622 2691
E-mail: info@consulab-bqsos.com

UNIPARTS S.A. DE C.V.
Tel: (0)55 5281 4718
Fax: (0)55 5281 4722
E-mail: uniparts@uniparts.com.mx

Netherlands

ENZO LIFE SCIENCES BVBA
Tel: +31/0 76 542 51 84
Fax: +31/0 76 542 52 61
E-mail: info-nl@enzolifesciences.com

New Zealand

SAPPHIRE BIOSCIENCE Pty. Ltd.
Tel: +61 (0) 2 9698 2022
Fax: +61 (0) 2 9698 1022
E-mail: sales@sapphirebioscience.com

UNITED BIORESEARCH PRODUCTS
Tel: +61 (0) 2 9651 3736
Fax: +61 (0) 2 9651 4247
E-mail: kirrily@unitedbioresearch.com.au

Norway

AH DIAGNOSTICS AS
Tel: (0)23 23 32 60
Fax: (0)23 23 32 70
E-mail: ahdiag@ahdiag.no

Pakistan

THE WORLDWIDE SCIENTIFIC
Tel: (0)42 755 2355
Fax: (0)42 755 3255
E-mail: wws@brain.net.pk

Poland

BIOMIBO
Tel: (0)22 872 07 97
Fax: (0)22 872 07 97
E-mail: biomibo@biomibo.com.pl

Portugal

BAPTISTA MARQUES, LDA
Tel: +351 (21) 722 06 60
Fax: +351 (21) 722 06 61
E-mail: geral@baptistamarques.pt

Romania

MEDIST SA
Tel: (0)21 411 5003
Fax: (0)21 410 5446
E-mail: office@medist.ro

Russia

CHIMMED Inc.
Tel: +7 095 728 4192
Fax: +7 095 742 8341
E-mail: bio@chimmed.ru

Singapore

ITS SCIENCE AND MEDICAL PTE. Ltd.
Tel: (0)6273 0898
Fax: (0)6273 0810
E-mail: info@its-sciencemedical.com

Slovakia

GENETICA s.r.o.
Tel: +42 (0) 2 7270 1055
Fax: +42 (0) 2 7270 1739
E-mail: genetika@genetica.cz

Slovenia

A-Ž LAB D.O.O.
Tel: +386 (0)1 433 63 22 /
+386 (0)1 230 18 84
Fax: +386 (0)1 230 19 85
E-mail: az.consulting@siol.net

South Africa

BIOCOM BIOTECH
Tel: +27 12 654 4614
Fax: +27 76 374 2093
E-mail: info@biocombiotech.co.za

Spain

GRUPO TAPER SA
Tel: +34 916 596 520
Fax: +34 916 610 084
E-mail: bioinvestigacion@grupotaper.com

Sweden

Immunkemi F&D AB
Tel: (0)8 583 615 00
Fax: (0)8 583 615 01
E-mail: sales@immunkemi.se

Switzerland

ENZO LIFE SCIENCES AG
Tel: +41 (0) 61 926 8989
Fax: +41 (0) 61 926 8979
E-mail: info-ch@enzolifesciences.com

Syria

NEW-MED TECHNOLOGY
Tel: (0)11 88271717
Fax: (0)11 88271710
E-mail: new-med@mail.sy

Taiwan

HONG JING Co., Ltd.
Tel: (0)2 3233 8585
Fax: (0)2 3233 8686
E-mail: hongjing6668@yahoo.com.tw

Thailand

THEERA TRADING CO., Ltd.
Tel: (0)2 412 5672 / (0)2 418 1068
Fax: (0)2 412 3244
E-mail: vtheera@ksc.th.com

ITS THAILAND CO., Ltd.
Tel: (0)2 308 0611
Fax: (0)2 308 0612
E-mail: info@its-thailand.com

Turkey

TOKRA MEDICAL Ltd.
Tel: (0)312 395 60 09
Fax: (0)312 395 39 61
E-mail: tokra@tokra.com.tr

Ukraine

CHIMMED Inc.
Tel: +7 095 728 4192
Fax: +7 095 742 8341
E-mail: bio@chimmed.ru

United Kingdom

ENZO LIFE SCIENCES (UK) Ltd.
Tel: 0845 601 1488 (UK customers)
Tel: +44 (0) 1392 825900 (from overseas)
Fax: +44 (0) 1392 825910
E-mail: info-uk@enzolifesciences.com

Uruguay

LAB SCIENTIFIC, INC. (USA)
Tel: (305) 716-9922
Fax: (305) 716-9923
E-mail: rshlesinger@labscent.com

USA

ENZO LIFE SCIENCES INTERNATIONAL, INC.
Tel: (610) 941-0430
Toll Free: 1-800-942-0430
Fax: (610) 941-9252
E-mail: info-usa@enzolifesciences.com

Venezuela

LAB SCIENTIFIC, INC. (USA)
Tel: (305) 716-9922
Fax: (305) 716-9923
E-mail: rshlesinger@labscent.com

Vietnam

ITS VIETNAM
Tel: (0)8 9255 232
Fax: (0)8 9255 233
E-mail: gmrs@its-vn.com

North/South America

ENZO LIFE SCIENCES INTERNATIONAL, INC.
5120 Butler Pike
Plymouth Meeting, PA 19462-1202 / USA
Tel. 1-800-942-0430 / (610) 941-0430
Fax (610) 941-9252
info-usa@enzolifesciences.com

Switzerland & Rest of Europe

ENZO LIFE SCIENCES AG
Industriestrasse 17, Postfach
CH-4415 Lausen / Switzerland
Tel. + 41/0 61 926 89 89
Fax + 41/0 61 926 89 79
info-ch@enzolifesciences.com

Benelux

ENZO LIFE SCIENCES BVBA
Melkerijweg 3
BE-2240 Zandhoven / Belgium
Tel. +32/0 3 466 04 20
Fax +32/0 3 466 04 29
info-be@enzolifesciences.com

France

ENZO LIFE SCIENCES FRANCE
c/o Covalab s.a.s
13, avenue Albert Einstein,
69100 Villeurbanne / France
Tel. +33/0 472 440 655
Fax +33/0 437 484 239
info-fr@enzolifesciences.com

Germany

ENZO LIFE SCIENCES GmbH
Marie-Curie-Strasse 8
DE-79539 Lörrach / Germany
Tel. +49/0 7621 5500 526
Fax +49/0 7621 5500 527
info-de@enzolifesciences.com

UK & Ireland

ENZO LIFE SCIENCES (UK) LTD.
Palatine House
Matford Court
Exeter EX2 8NL / UK
Tel. 0845 601 1488 (UK customers)
Tel. +44/0 1392 825900 (overseas)
Fax +44/0 1392 825910
info-uk@enzolifesciences.com

For local distributors see inside

