



## PERSONAL CARE



### Anti-Aging

Collagenases & Gelatinases

Proteases

Cell Proliferation

Oxidative Stress

### Irritation & Corrosion

### Allergy & Inflammation

### Genotoxicity

### Carcinogenesis

### Endocrine Disruption

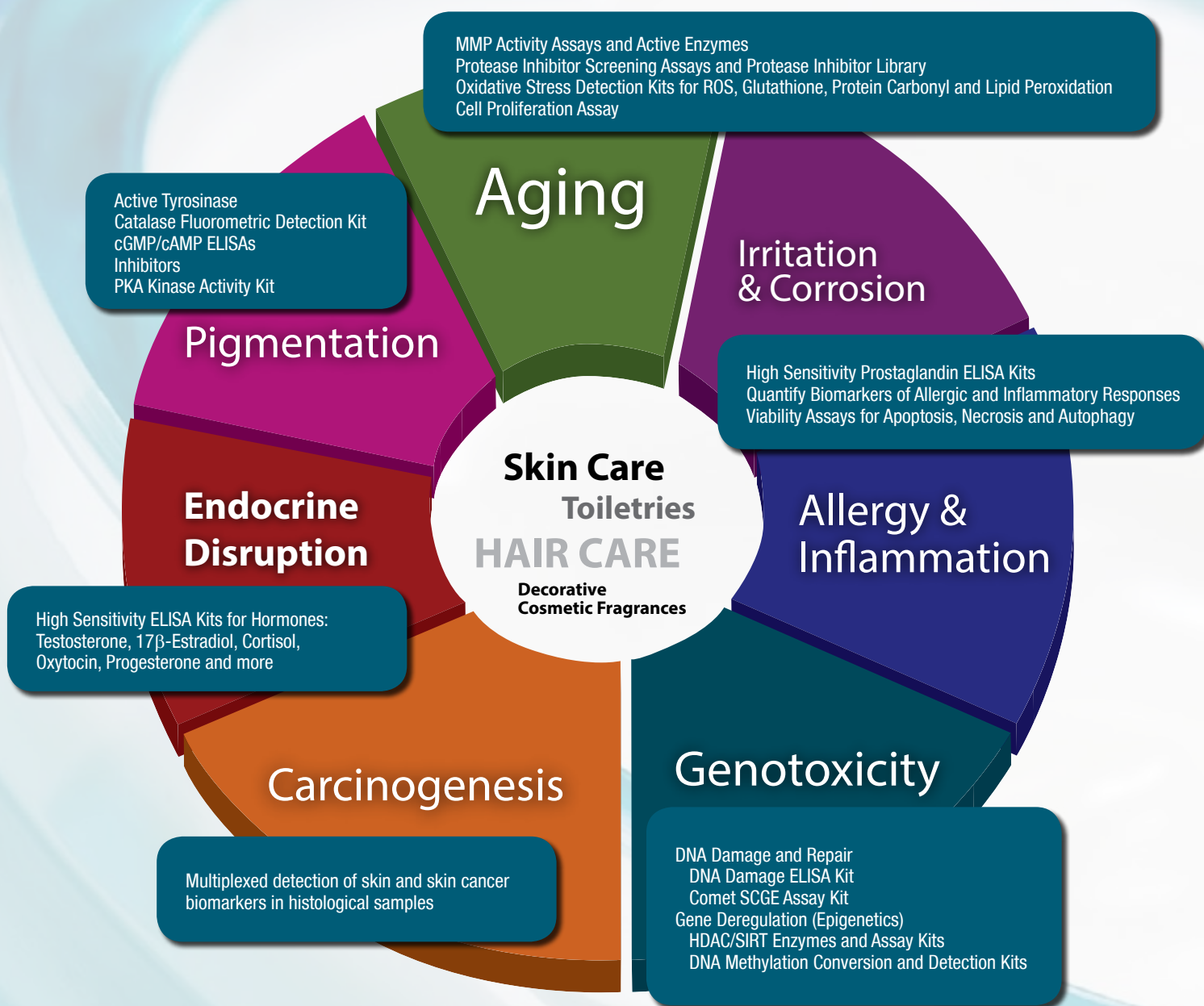
### Pigmentation

### Safety & Toxicity Testing

### Natural Products & Compound Screening

# INTEGRATED SOLUTIONS FOR PERSONAL CARE DEVELOPMENT

The skin functions as the primary line of defense against extrinsic stress such as UV-induced photo-damage, insults from microbial infections and physical deterioration resulting from aging and environmental exposure. As such, cosmetic industry scientists have utilized both basic and clinical research methods to develop effective ingredients for health and beauty products. The Enzo Life Sciences product portfolio provides assays and reagents for all phases of personal care product development to facilitate rapid discovery and enhanced testing for safety, toxicity and efficacy.



# CONTENTS

## Anti-Aging

### COLLAGENASES & GELATINASES

|  |   |
|--|---|
| Matrix Metalloproteinase (MMP) Inhibitor Profiling Kit | 4 |
| Matrix Metalloproteinase (MMP) Assay Kits              | 4 |
| Substrates   | 5 |
| Active Matrix Metalloproteinases (MMPs) and Proteins   | 6 |
| Inhibitors & Antibodies                                | 7 |

### PROTEASES

|   |   |
|---|---|
| Neutrophil Elastase Colorimetric Drug Discovery Kit | 8 |
| Protease Assay Kits & Substrates                    | 8 |
| Protease-Specific Antibodies                        | 8 |
| Active Proteases & Related Proteins                 | 9 |
| Protease Modulators                                 | 9 |

### CELL PROLIFERATION

|  |    |
|--|----|
| Cell Counting Kit                            | 10 |
| Proliferation-Related ELISA Kits             | 10 |
| Cell Proliferation Inhibitors                | 10 |
| NUCLEAR-ID® Cell Cycle Analysis Kits         | 10 |
| Growth Factors, Antibodies & Active Proteins | 11 |

### OXIDATIVE STRESS

|   |    |
|---|----|
| ROS Kits, Probes & Inhibitors                                 | 12 |
| Glutathione Detection Kits, Antibodies, Peptides & Modulators | 13 |
| Protein Carbonyl ELISA Kit                                    | 14 |
| Lipid Peroxidation Detection Kits & Modulators                | 14 |
| Cell Detoxification & Senescence                              | 15 |
| Proteostasis  | 15 |

## Irritation & Corrosion

|  |       |
|--|-------|
| Prostaglandin & Eicosanoid ELISA Kits                | 16    |
| Prostaglandin & Eicosanoid Antibodies & Biochemicals | 16-17 |
| Apoptosis & Necrosis Assay Kits                      | 17    |

## Allergy & Inflammation

|                                       |    |
|---------------------------------------|----|
| Pro-Inflammatory Cytokines ELISA Kits | 18 |
| Antibodies                            | 18 |

## Genotoxicity

|                                  |    |
|----------------------------------|----|
| DNA Damage ELISA Kit             | 19 |
| Comet SCGE Assay Kit             | 19 |
| Gene Dereglulation (Epigenetics) | 20 |

## Carcinogenesis

|                                     |       |
|-------------------------------------|-------|
| Immunohistochemistry (IHC) Reagents | 21    |
| IHC Antibodies                      | 22-23 |

## Endocrine Disruption

|                        |       |
|------------------------|-------|
| Hormone ELISA Kits     | 24-25 |
| Hormone Antibodies     | 26-27 |
| Biochemical Modulators | 28    |

## Pigmentation

|                            |    |
|----------------------------|----|
| cAMP ELISA Kits            | 29 |
| Inhibitors of Pigmentation | 29 |

## Safety & Toxicity Testing

|                     |    |
|---------------------|----|
| Cytotoxicity Assays | 30 |
|---------------------|----|

## Natural Products & Compound Screening

|  |    |
|--|----|
| Natural Product Library                    | 31 |
| Other Compound Libraries & Small Molecules | 31 |

## COLLAGENASES & GELATINASES

Collagenases (Matrix Metalloproteinase-1, MMP-8, and MMP-13) and gelatinases (MMP-2 and MMP-9) break down skin collagens, contributing to aging and wrinkling. Enzo offers high sensitivity colorimetric and fluorometric activity assays for screening collagenase and gelatinase inhibitors. Each assay is built using high purity active enzymes, and includes substrate, 96-well plate, and protocol.

### Optimized Detection of MMP Inhibitors

#### Matrix Metalloproteinase (MMP) Inhibitor Profiling Kit, Fluorometric RED (BML-AK308)

Inhibition of MMPs has been a target for inflammation, wound healing, skin aging, skin tone discoloration, rosacea, acne, and hair growth modulation. Our OMNIMMP® RED substrate offers key advantages over other substrates used for screening MMP inhibitors.

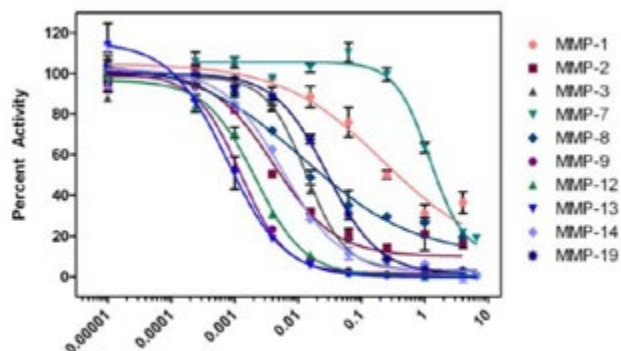
- Emission spectrum avoids the interference at lower wavelengths often exhibited by certain screening compounds
- Ultra-strong fluorescence allows for substrate concentrations much lower than the  $K_m$ , which is desirable for inhibitor screening assays
- Superior solubility characteristics compared with other MMP substrates
- Convenient real-time kinetics of cleavage is easily determined
- Includes active recombinant enzyme, substrate, and assay buffer

### Screen For Inhibitors of Specific MMPs

#### Matrix Metalloproteinase-1 (MMP-1) Fluorometric Drug Discovery Kit, RED (BML-AK301)

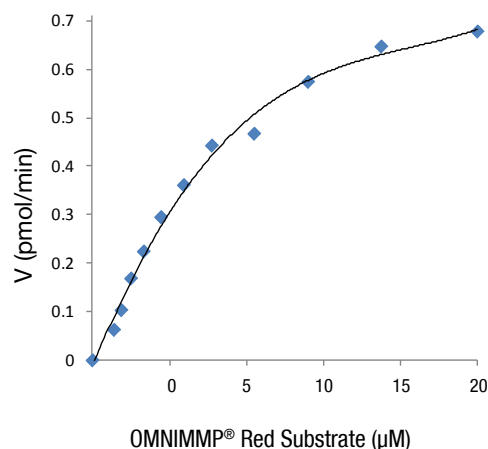
- Flexible platforms available for colorimetric and fluorometric detection
- Sensitive detection for specific broad range of MMPs
- Includes high quality substrate, enzyme, and prototypic control inhibitor

### Inhibitor Dose-response Curve



NNGH was pre-incubated with MMP enzyme before reactions were initiated by the addition of substrate.

### $K_{cat}/K_m$ Determination for MMP-1



| MMP Assay Kits   |           |            |
|--|-----------|------------|
| Product Name   | Product # | Size       |
| Matrix Metalloproteinase (MMP) Inhibitor Profiling Kit, Fluorometric Red | BML-AK308 | 1x96 Wells |
| MMP-1 Colorimetric Drug Discovery Kit                                    | BML-AK404 | 1x96 Wells |
| MMP-1 Fluorometric Drug Discovery Kit                                    | BML-AK405 | 1x96 Wells |
| MMP-1 Fluorometric Drug Discovery Kit, RED                               | BML-AK301 | 1x96 Wells |
| MMP-2 Colorimetric Drug Discovery Kit                                    | BML-AK408 | 1x96 Wells |

| MMP Assay Kits                                |           |            |
|---|-----------|------------|
| Product Name                                  | Product # | Size       |
| MMP-2 Fluorometric Drug Discovery Kit         | BML-AK409 | 1x96 Wells |
| MMP-2 Fluorometric Drug Discovery Kit, RED    | BML-AK302 | 1x96 Wells |
| MMP-3 Colorimetric Drug Discovery Kit         | BML-AK400 | 1x96 Wells |
| MMP-3 Fluorometric Drug Discovery Kit         | BML-AK401 | 1x96 Wells |
| MMP-3 Fluorometric Drug Discovery Kit, GREEN  | BML-AK303 | 1x96 Wells |
| MMP-3 Fluorometric Drug Discovery Kit, RED    | BML-AK311 | 1x96 Wells |
| MMP-7 Colorimetric Drug Discovery Kit         | BML-AK406 | 1x96 Wells |
| MMP-7 Fluorometric Drug Discovery Kit         | BML-AK407 | 1x96 Wells |
| MMP-7 Fluorometric Drug Discovery Kit, RED    | BML-AK304 | 1x96 Wells |
| MMP-8 Colorimetric Drug Discovery Kit         | BML-AK414 | 1x96 Wells |
| MMP-8 Fluorometric Drug Discovery Kit         | BML-AK415 | 1x96 Wells |
| MMP-8 Fluorometric Drug Discovery Kit, RED    | BML-AK305 | 1x96 Wells |
| MMP-9 Colorimetric Drug Discovery Kit         | BML-AK410 | 1x96 Wells |
| MMP-9 Fluorometric Drug Discovery Kit         | BML-AK411 | 1x96 Wells |
| MMP-9 Fluorometric Drug Discovery Kit, RED    | BML-AK306 | 1x96 Wells |
| MMP-10 Fluorometric Drug Discovery Kit        | BML-AK419 | 1x96 Wells |
| MMP-12 Colorimetric Drug Discovery Kit        | BML-AK402 | 1x96 Wells |
| MMP-12 Fluorometric Drug Discovery Kit        | BML-AK403 | 1x96 Wells |
| MMP-12 Fluorometric Drug Discovery Kit, GREEN | BML-AK312 | 1x96 Wells |
| MMP-13 Colorimetric Drug Discovery Kit        | BML-AK412 | 1x96 Wells |
| MMP-13 Fluorometric Drug Discovery Kit        | BML-AK413 | 1x96 Wells |
| MMP-14 Colorimetric Drug Discovery Kit        | BML-AK416 | 1x96 Wells |
| MMP-14 Fluorometric Drug Discovery Kit        | BML-AK417 | 1x96 Wells |
| MMP-19 Fluorometric Drug Discovery Kit, RED   | BML-AK307 | 1x96 Wells |
| MMP-20 Colorimetric Drug Discovery Kit        | BML-AK422 | 1x96 Wells |

| Substrates  |             |        |
|---|-------------|--------|
| Product Name  | Product #   | Size   |
| Cathepsin D & E Substrate (fluorogenic)                                     | BML-P145    | 1 mg   |
| Dnp-Pro-Leu-Gly-Met-Trp-Ser-Arg-OH  | ALX-260-120 | 5 mg   |
| Mca-Lys-Pro-Leu-Gly-Leu-Dpa-Ala-Arg-NH <sub>2</sub> (fluorogenic substrate) | BML-P276    | 1 mg   |
| Mca-Pro-Leu-Ala-Nva-Dap(Dnp)-Ala-Arg-NH <sub>2</sub>                        | ALX-260-123 | 1 mg   |
| Mca-RPPGFSAFK(Dnp)  | BML-P227    | 1 mg   |
| MMP Substrate (chromogenic)   | BML-P125    | 5 mg   |
| MMP Substrate (fluorogenic)   | BML-P128    | 1 mg   |
| MMP Substrate $\alpha$ (fluorogenic)  | BML-P131    | 1 mg   |
| MMP-3 Fluorogenic Substrate   | BML-P278    | 0.1 mg |
| OMNIMMP® Fluorogenic Control  | BML-P127    | 1 mg   |
| OMNIMMP® Fluorogenic Substrate  | BML-P126    | 1 mg   |
| OMNIMMP® RED Fluorogenic Substrate  | BML-P277    | 0.1 mg |
| TACE Substrate (fluorogenic)  | BML-P132    | 0.5 mg |

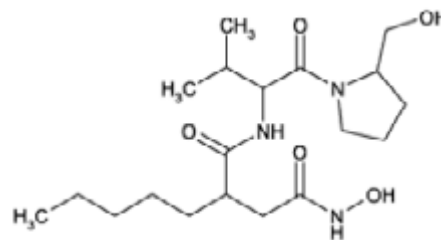
## ***Evaluation of extracellular matrix proteins and metalloproteinases***

Enzo provides a comprehensive portfolio of products for MMP research including active enzymes, antibodies, and high purity MMP inhibitors to suit your research needs.

| Active MMPs and Proteins   |             |            |
|--|-------------|------------|
| Product Name   | Product #   | Size       |
| Matrix Metalloproteinase (MMP) multipack-1                           | BML-AK013   | 10 µg each |
| Matrix Metalloproteinase (MMP) multipack-2                           | BML-AK014   | 10 µg each |
| MMP-1 (catalytic domain) (human), (recombinant)                      | BML-SE180   | 10 µg      |
| MMP-1 proenzyme (human fibroblasts)                                  | ALX-200-418 | 5 µg       |
| MMP-2 (catalytic domain) (human), (recombinant)                      | BML-SE237   | 10 µg      |
| MMP-2 (human fibroblasts), (purified)                                | BML-SE503   | 5 µg       |
| MMP-2 (mouse fibroblasts), (purified)                                | BML-SE505   | 5 µg       |
| MMP-2 proenzyme (human fibroblasts)                                  | ALX-200-419 | 5 µg       |
| MMP-2 proenzyme (human), (recombinant)                               | ALX-201-095 | 10 µg      |
| MMP-2/TIMP-2 proenzyme complex (human fibroblasts)                   | ALX-200-420 | 5 µg       |
| MMP-3 (catalytic domain) (human), (recombinant)                      | ALX-201-042 | 5 µg       |
| MMP-7 (catalytic domain) (human), (recombinant)                      | BML-SE181   | 10 µg      |
| MMP-7 (human), immunoblotting standard                               | BML-SW119   | 300 µL     |
| MMP-8 (catalytic domain) (human), (recombinant)                      | BML-SE255   | 10 µg      |
| MMP-8 (human neutrophils), (purified)                                | BML-SE492   | 5 µg       |
| MMP-8 proenzyme (human neutrophils)                                  | ALX-200-421 | 5 µg       |
| MMP-9 (catalytic domain) (human), (recombinant, <i>E. coli</i> )     | BML-SE360   | 10 µg      |
| MMP-9 (human neutrophils), (purified)                                | BML-SE504   | 5 µg       |
| MMP-9 (human neutrophils), dimer                                     | ALX-200-423 | 5 µg       |
| MMP-9 proenzyme (human neutrophils), monomer                         | ALX-200-422 | 5 µg       |
| MMP-9 proenzyme (human), monomer                                     | ALX-200-430 | 10 µg      |
| MMP-10 (catalytic domain) (human), (recombinant)                     | BML-SE329   | 10 µg      |
| MMP-11 (catalytic domain) (human), (recombinant)                     | BML-SE282   | 10 µg      |
| MMP-12 (catalytic domain), (human) (recombinant)                     | BML-SE138   | 10 µg      |
| MMP-13 (catalytic domain) (human), (recombinant)                     | BML-SE246   | 10 µg      |
| MMP-13 (human), (recombinant)  | BML-SE493   | 10 µg      |
| MMP-14 (catalytic domain) (human), (recombinant)                     | BML-SE259   | 10 µg      |
| MMP-14 (catalytic domain) (human), (recombinant)                     | ALX-201-098 | 10 µg      |
| MMP-14 prodomain (catalytic domain) (human), (recombinant) (His-tag) | ALX-201-099 | 10 µg      |
| MMP-14 proenzyme (soluble) (human), (recombinant) (His-tag)          | ALX-201-100 | 10 µg      |
| MMP-19 (catalytic domain) (human), (recombinant)                     | BML-SE561   | 10 µg      |
| MMP-20 (catalytic domain) (human), (recombinant)                     | BML-SE540   | 10 µg      |
| MMP-24 (catalytic domain) (human), (recombinant) (His-tag)           | ALX-201-105 | 10 µg      |

| MMP Inhibitors                      |             |             |
|-------------------------------------|-------------|-------------|
| Product Name                        | Product #   | Size        |
| Actinonin                           | ALX-260-128 | 5 mg        |
| CL-82198                            | BML-EI302   | 5 mg, 25 mg |
| CTT                                 | BML-PI136   | 1 mg, 5 mg  |
| Doxycycline . hyclate               | ALX-380-273 | 1 g, 5 g    |
| GM-6001                             | BML-EI300   | 1 mg, 5 mg  |
| Isobavachalcone                     | ALX-350-145 | 1 mg        |
| MMP-3 inhibitor                     | ALX-260-165 | 1 mg, 5 mg  |
| NNGH                                | BML-PI115   | 5 mg, 25 mg |
| PF-356231                           | BML-PI155   | 1 mg, 5 mg  |
| SB-3CT                              | BML-EI325   | 1 mg, 5 mg  |
| STT Negative Control for Gelatinase | BML-PI137   | 1 mg, 5 mg  |
| TAPI-0                              | BML-PI133   | 1 mg        |
| TAPI-1                              | BML-PI134   | 1 mg        |
| TAPI-2                              | BML-PI135   | 1 mg        |
| Z-Pro-Leu-Gly-NHOH                  | BML-PI139   | 100 mg      |

### Actinonin (ALX-260-128)

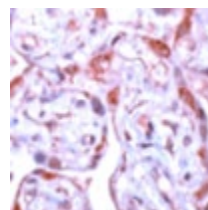


- Known inhibitor of select MMPs (-1, -2, -3, -7, -8, -9, -10, -12, -13)
- High purity, ≥98%
- Widely cited for various applications involving its inhibitory and antibiotic properties

| MMP-specific Antibodies            |              |        |
|------------------------------------|--------------|--------|
| Product Name                       | Product #    | Size   |
| MMP-1, pAb                         | ADI-905-472  | 1 mL   |
| MMP-1, pAb                         | ALX-210-876  | 100 µg |
| MMP-2 (CT), pAb                    | BML-SA616    | 50 µg  |
| MMP-2 (hinge region), pAb          | BML-SA679    | 100 µg |
| MMP-2, pAb                         | ALX-210-753  | 500 µL |
| MMP-3, pAb                         | BML-SA104    | 100 µg |
| MMP-7 (CT), pAb                    | BML-SA625    | 50 µg  |
| MMP-7, mAb (2D3)                   | ADI-905-1013 | 100 µg |
| MMP-9 (CT), pAb                    | BML-SA620    | 500 µL |
| MMP-9 (hinge region), pAb          | BML-SA680    | 100 µg |
| MMP-9, pAb                         | ALX-210-756  | 100 µg |
| MMP-10 (human) (hinge region), pAb | BML-SA434    | 100 µg |
| MMP-12 (CT), pAb                   | BML-SA467    | 100 µL |
| MMP-12 (hinge region), pAb         | BML-SA669    | 100 µg |
| MMP-13 (hinge region), pAb         | BML-SA626    | 50 µg  |
| MMP-13 (human), mAb (M53)          | ALX-804-365  | 100 µL |
| MMP-13 (human), pAb                | ALX-210-833  | 100 µL |

### MMP-1, pAb (ADI-905-472)

- Recognizes human MMP-1
- Validated for IHC



Immunohistochemical analysis of human placenta stained with MMP-1, pAb.



## PROTEASES

### *Rapidly screen inhibitors of key enzymes regulating elasticity*

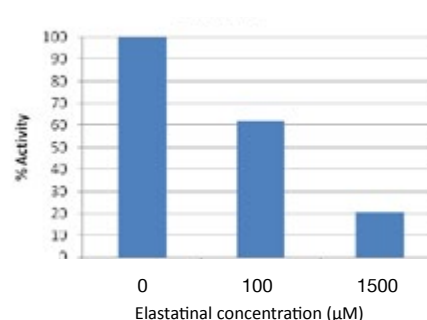
Cathepsins, Granzymes and Elastases are involved in regulatory mechanisms in human skin and act as biomarkers for decreased skin elasticity due to UV exposure in photoaged human skin.

#### **Neutrophil Elastase Colorimetric Drug Discovery Kit (BML-AK497)**

Neutrophil elastase is a serine protease found in polymorphonuclear neutrophils that functions in inflammation.

- Rapid, sensitive, and reliable results for the screening of neutrophil elastase inhibitors
- Kit includes elastatinal, a known elastase inhibitor, as a control
- Amenable to high-throughput screening

### **Percent Inhibition of HNE by Elastatinal**



Using the elastase inhibitor, elastatinal, the % activity remaining can be calculated.

| Protease Assay Kits & Substrates                    |             |              |
|---|-------------|--------------|
| Product Name  | Product #   | Size         |
| Cathepsin D & E Substrate (fluorogenic)             | BML-P145    | 1 mg         |
| Cathepsin K Drug Discovery Kit                      | BML-AK430   | 1x96 Wells   |
| Cathepsin S Drug Discovery Kit                      | BML-AK431   | 1x96 Wells   |
| CV-Cathepsin B Detection Kit                        | BML-AK125   | 1x96 Wells   |
| CV-Cathepsin K Detection Kit                        | BML-AK126   | 1x96 Wells   |
| CV-Cathepsin L Detection Kit                        | BML-AK127   | 1x96 Wells   |
| Granzyme B Assay Kit for Drug Discovery             | BML-AK711   | 1x96 Wells   |
| Neutrophil Elastase Colorimetric Drug Discovery Kit | BML-AK497   | 1x96 Wells   |
| OMNICATHEPSIN® Fluorogenic Substrate                | BML-P139    | 10 mg, 50 mg |
| PMN-Elastase (Human) ELISA Kit                      | ALX-850-265 | 1x96 Wells   |

| Protease-Specific Antibodies                       |                |              |
|--|----------------|--------------|
| Cathepsin B (human), pAb                           | BML-SA361      | 100 μL       |
| Cathepsin D (human), pAb                           | BML-SA398      | 100 μL       |
| Cathepsin G (human), pAb                           | BML-SA354      | 100 μL       |
| Cathepsin H (human), pAb                           | BML-SA355      | 100 μL       |
| Cathepsin L (human), pAb                           | BML-SA362      | 100 μL       |
| Cathepsin L/procathepsin L (human), mAb (CLP 1/36) | ALX-804-292    | 200 μg       |
| Cathepsin L/procathepsin L, mAb (CPLH 33/2)        | ALX-804-291    | 100 μg       |
| Cathepsin L/procathepsin L, mAb (CPLH 3G10)        | ALX-804-293    | 200 μg       |
| Collagen Type I, mAb (5D8)                         | BPD-CSI-008-01 | 200 μg, 1 mg |
| Collagen Type III, mAb (1E7-D7/Col3)               | BPD-CSI-007-01 | 200 μg       |
| Collagen Type XII, mAb (378D5)                     | ALX-803-309    | 100 μL       |
| Granzyme B (human), mAb (B18.1)                    | ALX-804-121    | 100 μg       |
| Granzyme B (human), mAb (B18.1) (FITC conjugate)   | ALX-804-121F   | 100 Tests    |
| Granzyme B (human), mAb (GrB7)                     | ALX-804-198    | 50 μg        |
| Neutrophil elastase (human), (purified)            | BML-SE284      | 100 μL       |

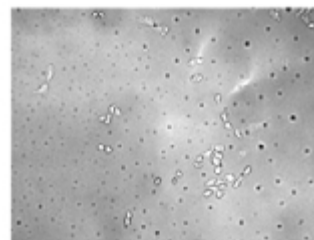


| Active Proteases and Related Proteins           |             |               |
|---|-------------|---------------|
| Product Name                                    | Product #   | Size          |
| Cathepsin B (human liver), (purified)           | BML-SE198   | 25 µg         |
| Cathepsin D (human liver), (purified)           | BML-SE199   | 25 µg         |
| Cathepsin D (human), (recombinant) (His-tag)    | ADI-908-241 | 10 µg         |
| Cathepsin F (human), (recombinant)              | BML-SE568   | 5 µg          |
| Cathepsin G (human neutrophils), (purified)     | BML-SE283   | 100 µg        |
| Cathepsin H (human liver), (purified)           | BML-SE200   | 25 µg         |
| Cathepsin K (human), (recombinant)              | BML-SE553   | 10 µg         |
| Cathepsin L (human liver), (purified)           | BML-SE201   | 25 µg         |
| Cathepsin S (human), (recombinant)              | BML-SE453   | 10 µg         |
| Collagen I, rat tail                            | ALX-522-435 | 20 mg, 100 mg |
| Collagen I, rat tail (thin plate coating)       | ALX-522-440 | 50 mL         |
| Elafin (human), (recombinant)                   | ALX-201-240 | 100 µg        |
| Granzyme A (human)                              | ALX-200-605 | 10 µg         |
| Granzyme A (human), (recombinant)               | ALX-201-118 | 10 µg         |
| Granzyme B (human lymphocytes)                  | ALX-200-602 | 10 µg         |
| Granzyme B (human), (recombinant)               | BML-SE238   | 5000 U        |
| Granzyme K (human), (recombinant)               | ALX-201-117 | 10 µg         |
| Neutrophil elastase (human), (purified)         | BML-SE284   | 100 µg        |
| Procathepsin F (human), (recombinant)           | BML-SE541   | 5 µg          |
| Procathepsin K (human), (recombinant)           | BML-SE367   | 10 µg         |
| Procathepsin K (human), (recombinant)           | ALX-201-239 | 10 µg         |
| Procathepsin V (human), (recombinant) (His-tag) | BML-SE554   | 10 µg         |
| SLPI (human), (recombinant) (His-tag)           | ADI-908-304 | 10 µg         |

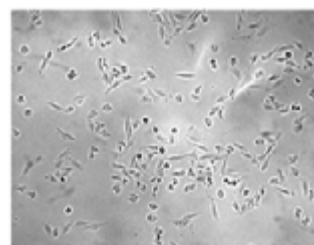
### Collagen I, Rat Tail (ALX-522-435)

- High Purity, ≥90% (SDS-PAGE)
- Sterile format for cell culture applications
- Also available as a ready-to-use reagent for thin plate coating (ALX-522-440)

#### Untreated



#### Treated



Collagen I was tested in a cell proliferation assay. Increased attachment of cells on collagen-coated coverslips was observed following culture.

| Protease Modulators     |           |             |
|-------------------------|-----------|-------------|
| Product Name            | Product # | Size        |
| 3,4 Dichloroisocoumarin | BML-PI110 | 10 mg       |
| Azelastine . HCl        | BML-NH108 | 100 mg      |
| Elastatinal             | BML-PI103 | 5 mg, 25 mg |

## Are you screening for protease inhibitors?

Our SCREEN-WELL® Protease Inhibitor Library includes 53 known protease inhibitors on a single 96-well plate.

## CELL PROLIFERATION

### Convenient Assays, Sensitive Results

#### Cell Counting Kit-8 (ALX-850-039)

A high-sensitivity colorimetric assay for the determination of the number of viable cells, using WST-8, in cell proliferation and cytotoxicity assays. This critical assay can be used to screen modulators of cell proliferation that affect the appearance of wrinkles and other signs of aging.

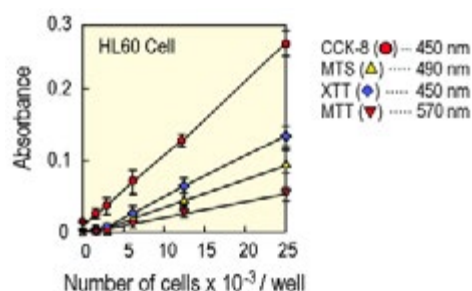
- More sensitive than MTT, XTT, MTS or WST-1
- Ready-to-use solution can be added directly to the cell media
- No organic solvents or isotopes required
- No harvesting, washing, or solubilization steps
- Amenable to high-throughput screening

| Cell Proliferation Assay |             |                        |
|--------------------------|-------------|------------------------|
| Product Name             | Product #   | Size                   |
| Cell Counting Kit-8      | ALX-850-039 | 500 Tests, 5x500 Tests |

| Proliferation Related ELISA Kits                           |              |            |
|--|--------------|------------|
| Product Name   | Product #    | Size       |
| Erk1/2 ELISA Kit   | ADI-900-152  | 1x96 Wells |
| [pThr <sup>202</sup> /Tyr <sup>204</sup> ]Erk1/2 ELISA Kit | ADI-900-098A | 1x96 Wells |
| IGF-1 (human), ELISA Kit                                   | ADI-900-150  | 1x96 Wells |
| TGF-β1 ELISA Kit   | ADI-900-155  | 1x96 Wells |

| Cell Proliferation Inhibitors |             |            |
|-------------------------------|-------------|------------|
| Product Name                  | Product #   | Size       |
| Tyrphostin AG 1024            | ALX-270-217 | 1 mg, 5 mg |
| Reveromycin A                 | ALX-380-216 | 0.25 mg    |
| Reveromycin D                 | ALX-380-219 | 0.25 mg    |

### High-sensitivity Cell Proliferation Assay



CTLL-2 cells were incubated with various concentrations of IL-2 for 72 hours. CCK-8 solution was added to each well and the absorbance at 450 nm was measured. IL-2 exposure resulted in increased absorbance, which correlates to an increase in cell proliferation.

### The Brightest, Most Sensitive Cell Cycle Analysis Dye for Live Cells

#### NUCLEAR-ID® Red Cell Cycle Kit (ENZ-51008)

Our stable NUCLEAR-ID® Red DNA stain crosses the membrane of live cells and specifically targets dsDNA. The dye can be used for various applications including nucleated cell gating in flow cytometry, DNA ploidy and cell cycle analysis, and cytoplasm segmentation in high content screening (HCS).

- Brighter, photostable red dye yields improved sensitivity
- Highly cell-permeable dye minimizes cell density optimization
- Lower cost per test than comparable dyes
- Suitable for multiplexing with green fluorophores, including GFP and FITC
- Available as a complete kit with controls or as a stand-alone dye

| NUCLEAR-ID® Cell Cycle Analysis Kits            |           |                    |
|---|-----------|--------------------|
| Product Name                                    | Product # | Size               |
| NUCLEAR-ID® Red Cell Cycle Kit (GFP-CERTIFIED®) | ENZ-51008 | 1 Kit (100 assays) |
| NUCLEAR-ID® Green Cell Cycle Kit                | ENZ-51014 | 1 Kit (100 assays) |
| NUCLEAR-ID® Red DNA Stain                       | ENZ-52406 | 200 µL             |

### Brightest Dye for Cell Cycle Analysis



## Growth Factors

Growth factors are frequently used in cosmetics to reverse the signs and symptoms of aging skin mediated by environmental elements.

| Antibodies                                       |             |        |
|--|-------------|--------|
| Product Name                                     | Product #   | Size   |
| Fibroblast growth factor 2, mAb (3D9)            | ADI-905-832 | 100 µL |
| Fibroblast growth factor 2, mAb (2H11)           | ADI-905-833 | 100 µL |
| Fibroblast growth factor 1, mAb (4D2)            | ADI-905-883 | 100 µL |
| Insulin-like growth factor-1 receptor, mAb (1-2) | BML-SA369   | 100 µg |
| TGF-βRIII (human), mAb (Jack-1)                  | ALX-804-871 | 100 µg |

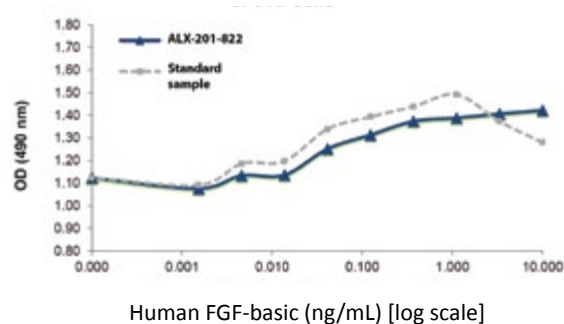
| Active Proteins  |             |               |
|--|-------------|---------------|
| Product Name   | Product #   | Size          |
| EGF (human), (recombinant)   | ADI-908-060 | 100 µg        |
| Epidermal Growth Factor (rat), (recombinant)                           | ALX-201-812 | 100 µg        |
| FGF-21 (mouse), (recombinant)  | ALX-201-401 | 10 µg         |
| FGF-21 (mouse), (recombinant) (His-tag)                                | ALX-201-409 | 10 µg         |
| FGF-basic 147 (human), (recombinant)                                   | ALX-201-822 | 10 µg, 50 µg  |
| FGF-basic 154 (human), (recombinant)                                   | ALX-201-823 | 10 µg, 50 µg  |
| FGF basic (mouse), (recombinant)                                       | ALX-201-824 | 10 µg, 50 µg  |
| IGF-1 (human), (recombinant)   | ADI-908-059 | 20 µg         |
| Insulin-like Growth Factor-1 Receptor (human), (recombinant) (GST-tag) | BML-SE232   | 20 µg         |
| TGF-β3 (mouse), (recombinant)  | ALX-201-817 | 10 µg, 100 µg |

### FGF-basic 147 (human) (ALX-201-822)

A critical factor in angiogenesis, wound healing and embryonic development

- Carrier-free
- High purity, ≥ 97%
- Activity tested in cell proliferation assay

### Human FGF-basic-induced Proliferation of 3T3 Cells



3T3 cells were cultured with 0 to 10 ng/mL human FGF-basic. Cell proliferation was measured after 44 hours and the linear portion of the curve was used to calculate the ED50. The ED50 for this lot of human FGF-basic was 0.042-0.063 ng/mL. The typical expected range is less than 1 ng/mL.

# ANTI-AGING

## Determination of Antioxidant Capacity Through Oxidative Stress Biomarkers

Oxidative stress in response to aging and environmental stimuli can lead to tissue damage. Analysis of reactive oxygen species, mitochondrial dysfunction, lipid peroxidation and glutathione levels can provide an indication of the effectiveness of cosmetic and cosmeceutical formulations. Enzo offers a selection of assay kits which enable detection of oxidative stress from the nucleic acid level up through a complete whole cell analysis.

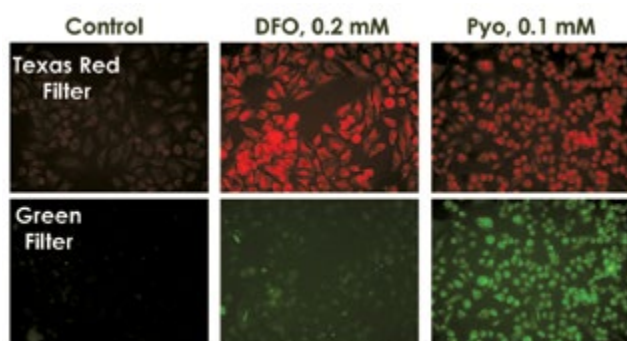
## REACTIVE OXYGEN SPECIES

### Multiplex, Real-time Analysis of Hypoxia and ROS in Live Cells

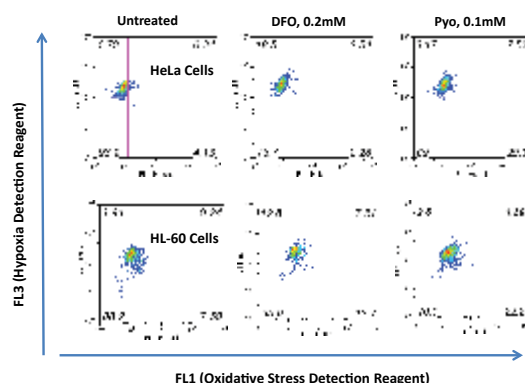
#### ROS-ID® Hypoxia/Oxidative Stress Detection Kit (ENZ-51042)

- Sensitive hypoxia dye fluoresces red when converted by nitroreductases
- Total ROS dye fluoresces green in presence of a variety of reactive oxygen species
- Non-toxic, cell-permeable dyes optimized to avoid artifactual fluorescence
- Suitable for microscopic or flow cytometric analysis of adherent or suspension cells
- Amenable to high-throughput screening










#### Simultaneously Monitor Hypoxic Status and ROS Formation



#### Quantitative Multiparametric Assay for Flow Cytometry



Bright red fluorescence of the hypoxia probe is observed following its conversion by cellular nitroreductases under hypoxic conditions such as those induced chemically by treatment with the hypoxia-mimetic desferrioxamine (DFO). The assay facilitates simultaneous quantification of hypoxic status (FL3/red) and presence of ROS (FL1/green) by flow cytometry.

| ROS Detection Kits                             |           |   |   |
|--|-----------|---|---|
| Product Name                                   | Product # | Application   | Size  |
| ROS-ID® Hypoxia/Oxidative Stress Detection Kit | ENZ-51042 |     | 500 Microscopy or 100 Flow Cytometry Assays |
| ROS-ID® ROS/RNS Detection Kit for Microscopy   | ENZ-51001 |    | 200 Assays                                  |
| ROS-ID® Total ROS/Superoxide Detection Kit     | ENZ-51010 |    | 200 Microscopy or 50 Flow Cytometry Assays  |
| ROS-ID® Total ROS Detection Kit                | ENZ-51011 |    | 200 Microscopy or 50 Flow Cytometry Assays  |

| Probes for ROS                    |             |           |
|-----------------------------------|-------------|-----------|
| Product Name                      | Product #   | Size      |
| 5(6)-CDCFDA (ultra pure)          | ENZ-52103   | 100 mg    |
| CFDA (ultra pure)                 | ENZ-52104   | 100 mg    |
| Dihydrorhodamine 123 (ultra pure) | ENZ-52302   | 10 mg     |
| hROS, (detection reagent)         | ADI-906-043 | 150 Tests |
| APF                               | ALX-620-075 | 1 mg      |
| HPF                               | ALX-620-074 | 1 mg      |

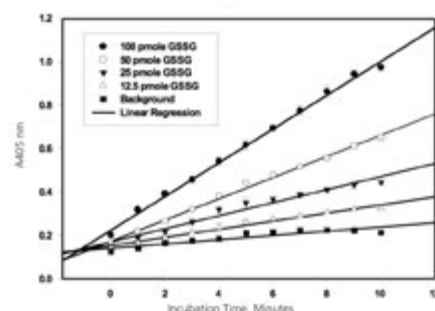
# GLUTATHIONE

## Glutathione (Total) Detection Kit (ADI-900-160)

Glutathione Reductase reduces oxidized glutathione (GSSG) to reduced glutathione (GSH). The sulfhydryl group of GSH reacts with DTNB (5,5'-dithiobis-2-nitrobenzoic acid, Ellman's reagent) to produce a yellow colored 5-thio-2-nitrobenzoic acid (TNB) that absorbs at 405 or 414 nm. The rate of TNB production is directly proportional to the concentration of glutathione in the sample.

- High-sensitivity enzymatic recycling reaction measures total, oxidized, or reduced glutathione
- Broad detection range: range (12.5 - 100 pmol/well)
- Easy-to-use sample handling protocols for a wide variety of matrices
- Convenient colorimetric 96-well plate format

## Measure Oxidized and/or Reduced Glutathione



Change in absorbance at 405 nm versus incubation time as a function of pmoles of GSSG/well.

## Glutathione Assay Kits & Probes

| Product Name                        | Product #   | Size       |
|-------------------------------------|-------------|------------|
| Glutathione (total), Detection Kit  | ADI-900-160 | 4x96 Wells |
| Glutathione Peroxidase Activity Kit | ADI-900-158 | 5x96 Wells |
| Glutathione Reductase Activity Kit  | ADI-900-159 | 5x96 Wells |
| Monobromobimane (ultra pure)        | ENZ-52501   | 25 mg      |

## Glutathione Antibodies

| Product Name                            | Product #   | Size   |
|---|-------------|--------|
| Glutathione, mAb (D8)                   | ADI-SPA-542 | 100 µg |
| Glutathione S-Transferase, mAb (8C1-G4) | ADI-905-618 | 100 µg |

## Glutathione Chemical Modulators

| Product Name                                   | Product #   | Size                 |
|--|-------------|----------------------|
| 5,8,11-Eicosatriynoic acid                     | BML-ET003   | 1 mg, 10 mg          |
| Butein   | ALX-350-246 | 10 mg                |
| Buthionine sulfoximine                         | BML-FR117   | 500 mg               |
| Cafestol                                       | ALX-350-220 | 50 mg                |
| Carmustine                                     | ALX-400-037 | 25 mg                |
| Curcumin (high purity)                         | ALX-350-028 | 10 mg, 50 mg, 250 mg |
| Ebselen  | ALX-270-097 | 5 mg, 25 mg          |
| Ethacrynic acid                                | BML-EI128   | 1 g                  |
| Kahweol  | ALX-350-223 | 10 mg                |
| Prostaglandin A2                               | BML-PG002   | 1 mg, 10 mg          |
| RSSR   | ALX-430-102 | 10 mg, 25 mg, 50 mg  |
| S-(N-(3-Phenylpropyl)thiocarbamoyl)-L-cysteine | BML-T114    | 50 mg                |
| Sedanolid                                      | ALX-350-229 | 100 mg               |

## PROTEIN CARBOXYLATION

***Protein carbonylation is a known irreversible result of oxidative damage that can lead to protein aggregation and dysfunction***

### **Protein Carbonyl ELISA Kit (ALX-850-312)**

This convenient ELISA kit enables quantitative determination of carbonylated protein levels in plasma, other body fluids, cell and tissue extracts.

- High-sensitivity
- Low background
- Amenable to high-throughput analysis
- Validated and widely cited for various sample matrices

## LIPID PEROXIDATION

Lipid peroxidation is a well-known example of oxidative damage, and lipid peroxides are prominent non-radical intermediates of lipid peroxidation. We offer a variety of assays and small molecules for screening the effects of anti-oxidants as potential inhibitors of this pathway.

### **ALDetect™ (MDA-specific) Lipid Peroxidation Assay Kit (BML-AK171)**

Designed to assay free malondialdehyde (MDA) or, after a hydrolysis step, total MDA (i.e., free and protein-bound Schiff base conjugates). Features include:

- Minimal interference from other lipid peroxidation products, such as 4-hydroxyalkenals
- Assess MDA levels from a variety of sample types including tissue and cell lysates
- Rapid results in less than 1.5 hours

| Lipid Peroxidation Assays                             |             |           |
|---|-------------|-----------|
| Product Name  | Product #   | Size      |
| ALDetect™ Lipid Peroxidation assay Kit                | BML-AK170   | 100 Tests |
| ALDetect™ (MDA-specific) Lipid Peroxidation assay Kit | BML-AK171   | 25 Tests  |
| OXI-TEK TBARS Assay Kit                               | ALX-850-287 | 160 Tests |

| Biochemical Modulators of Oxidative Stress |             |                   |
|--|-------------|-------------------|
| (-)-Epigallocatechin gallate               | ALX-270-263 | 10 mg, 50 mg      |
| (Z)-4-Hydroxytamoxifen                     | ALX-550-361 | 1 mg, 5 mg        |
| Auraptene                                  | ALX-350-361 | 5 mg, 25 mg       |
| Bakuchiol                                  | ALX-350-144 | 1 mg              |
| Caffeic acid phenylethyl ester             | ALX-270-244 | 10 mg, 50 mg      |
| Carazostatin                               | ALX-350-253 | 100 µg, 1 mg      |
| Carnosic acid                              | ALX-270-264 | 10 mg, 50 mg      |
| Celastrol                                  | ALX-350-332 | 5 mg, 25 mg       |
| Cinnamtannin B-1                           | ALX-350-365 | 5 mg              |
| Idebenone                                  | BML-FR114   | 20 mg, 100 mg     |
| Morin                                      | ALX-385-016 | 1 g               |
| Ochratoxin A                               | ALX-630-089 | 1 mg, 5 mg, 25 mg |
| Pyrrolostatin                              | ALX-350-252 | 100 µg, 1 mg      |
| Rosmarinic acid                            | ALX-270-253 | 10 mg, 50 mg      |
| Trolox®                                    | ALX-270-267 | 100 mg            |
| U-74389G                                   | BML-FR100   | 100 mg            |

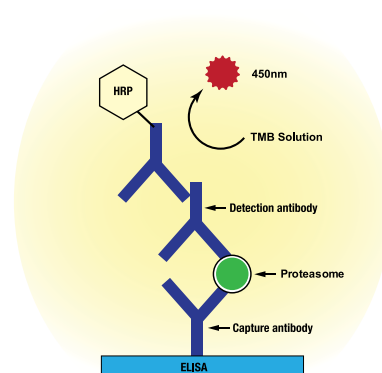
## CELL DETOXIFICATION

**Examine the link between cell detoxification and proteasome concentrations with the only commercially available kit for proteasome quantification**

### Proteasome ELISA Kit (BML-PW0575)

- Quantify 20S proteasome concentrations in biological samples (serum, plasma, cell lysates, tissue extracts)
- Investigate effects of inhibitors and activations on levels of circulating proteasome

### Quantify Circulating Levels of 20S Proteasome



Summary of Proteasome ELISA protocol

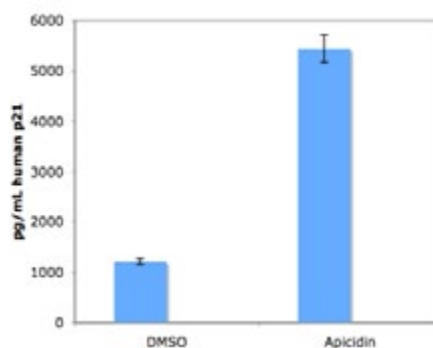
## CELL SENEESCENCE

### p21 ELISA Kit (ADI-900-161)

The p21 (human) ELISA kit is a colorimetric, immunometric enzyme immunoassay kit with results in 3 hours. Save time, money, and precious sample with fully quantitative results and increased sensitivity compared to Western blot analysis.

- Sensitive ELISA quantifies picogram levels compared to microgram levels in Western blot
- Excellent lot-to-lot reproducibility
- Easy-to-use liquid color-coded reagents reduce error

### Monitor Levels of p21, a Regulator of Senescence



Levels of p21 in control (DMSO- or Apicidin-treated HeLa cell lysates) as determined using p21 ELISA kit.

| Related Products                             |             |            |
|--|-------------|------------|
| Product Name                                 | Product #   | Size       |
| p21 ELISA Kit                                | ADI-900-161 | 1x96 Wells |
| Cellular Senescence Activity Assay           | ENZ-KIT129  | 120 Assays |
| Cellular Senescence Live Cell Analysis Assay | ENZ-KIT130  | 10 Assays  |

## PROTEOSTASIS

### AMP'D® HSP70 High Sensitivity ELISA Kit (ENZ-KIT101)

Ultra-sensitive AMP'D® HSP70 high sensitivity ELISA kit enabling the ability to use less sample and detect both baseline and upregulated levels of human, mouse and rat Hsp70 (Hsp72), a major chaperone, cancer biomarker, and key cell stress regulator.

- Ultra-sensitive measurement of Hsp70, detecting as little as 7 pg/mL
- Negligible reactivity from similar Hsp70 family members (Hsc70/Hsp73, Grp78, DnaK, or Hsp71)
- High-throughput format with results in 4.5 hours for up to 38 samples in duplicate
- Fully quantitative results that surpass semi-quantitative Western blot analysis

| Related Products                 |              |            |
|----------------------------------|--------------|------------|
| Product Name                     | Product #    | Size       |
| HSP70 ELISA Kit                  | ADI-EKS-700B | 1x96 Wells |
| HSP70 High Sensitivity ELISA Kit | ADI-EKS-715  | 1x96 Wells |



# IRRITATION & CORROSION

## PROSTAGLANDINS AND OTHER EICOSANOIDS

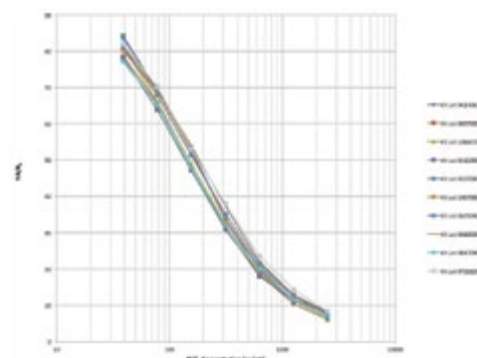
### *All of the sensitivity. None of the doubt.*

#### Highest Sensitivity, Widely Cited PGE<sub>2</sub> ELISA Kits

For over two decades, scientists around the world have used Enzo's ELISA kits to reliably detect PGE<sub>2</sub> in their samples. Over that time, we have continued to improve our kits to enable PGE<sub>2</sub> detection with better sensitivity, flexibility, and consistency.

- Ultrasensitive ELISAs to measure as little as 8.26 pg/mL PGE<sub>2</sub>
- Compatible with culture supernatants, serum, saliva, urine, whole blood and more
- Widely cited in peer reviewed literature for over two decades
- High-throughput assays in colorimetric, chemiluminescent, and fluorescent readouts
- Time-tested manufacturing provides consistent lot-to-lot results

### Reliable Manufacturing Ensures Reproducibility



Graph demonstrating the robust and reproducible nature of PGE<sub>2</sub> ELISA Kits showing standard curves from ten kit lots manufactured over five years.

| Prostaglandin & Eicosanoids ELISA Kits       |             |             |
|--|-------------|-------------|
| Product Name                                 | Product #   | Size        |
| 8-iso-PGF <sub>2α</sub> ELISA Kit            | ADI-900-010 | 1x96 Wells* |
| 11-dehydro-TXB <sub>2</sub> ELISA Kit        | ADI-900-092 | 1x96 Wells* |
| 12(S)-HETE ELISA Kit                         | ADI-900-050 | 1x96 Wells* |
| 15(S)-HETE ELISA Kit                         | ADI-900-051 | 1x96 Wells* |
| Cysteinyl leukotriene ELISA Kit              | ADI-900-070 | 1x96 Wells* |
| Direct 8-iso-PGF <sub>2α</sub> ELISA Kit     | ADI-900-091 | 1x96 Wells* |
| LTB <sub>4</sub> ELISA Kit                   | ADI-900-068 | 1x96 Wells* |
| PGE <sub>2</sub> CLIA Kit                    | ADI-910-001 | 1x96 Wells* |
| PGE <sub>2</sub> ELISA Kit                   | ADI-900-001 | 1x96 Wells* |
| PGE <sub>2</sub> FPIA Kit                    | ADI-920-001 | 100 Tests   |
| PGE <sub>2</sub> High Sensitivity ELISA Kit  | ADI-930-001 | 1x96 Wells* |
| PGF <sub>2α</sub> ELISA Kit                  | ADI-900-069 | 1x96 Wells* |
| PGF <sub>2α</sub> High Sensitivity ELISA Kit | ADI-930-069 | 1x96 Wells* |
| TXB <sub>2</sub> ELISA Kit                   | ADI-900-002 | 1x96 Wells* |

| Prostaglandin & Eicosanoids Antibodies |             |        |
|--|-------------|--------|
| Product Name                           | Product #   | Size   |
| 12(L)-HETE, pAb                        | ADI-905-053 | 100 µL |
| 8-iso-PGF <sub>2α</sub> , pAb          | ADI-905-015 | 100 µL |
| 8-iso-PGF <sub>2α</sub> , pAb          | ADI-905-016 | 100 µL |
| PGE <sub>2</sub> , mAb (BG8)           | ADI-905-628 | 100 µL |
| PGE <sub>2</sub> , pAb                 | ADI-905-013 | 100 µL |
| PGE <sub>2</sub> , pAb                 | ADI-905-025 | 1 mL   |
| TXB <sub>2</sub> , pAb                 | ADI-905-005 | 100 µL |

\*5x96 well pack size also available

| Related Biochemicals  |             |               |
|---|-------------|---------------|
| Product Name  | Product #   | Size          |
| (±)13-Azaprostanoic acid                                      | BML-PG036   | 1 mg, 10 mg   |
| 12(S)-Hydroperoxyeicosa-5Z,8Z,10E,14Z-tetraenoic acid (HPETE) | BML-HP012   | 50 µg         |
| Caffeic acid ethyl ester                                      | ALX-270-480 | 50 mg, 250 mg |
| Carbethoxyhexyl imidazole                                     | BML-EI120   | 10 mg, 50 mg  |
| Carboxyheptyl imidazole                                       | BML-EI119   | 10 mg, 50 mg  |
| Cysteinyl leukotriene mixture                                 | BML-SM105   | 10 µg each    |
| Disodium cromoglycate   | BML-EI121   | 1 g, 5 g      |
| Eicosapentaenoic acid   | BML-FA001   | 100 mg, 1 g   |
| Furegrelate sodium  | ALX-270-120 | 10 mg, 50 mg  |
| Ketoconazole  | BML-EI107   | 50 mg, 500 mg |
| L-655,240   | BML-RA117   | 1 mg, 5 mg    |
| Leukotriene B4 & ω oxidation metabolites mixture              | BML-SM115   | 10 µg each    |
| Ozagrel   | BML-EI378   | 10 mg         |
| Picotamide  | BML-RA104   | 10 mg, 50 mg  |
| Seratrodast   | BML-RA119   | 10 mg, 50 mg  |
| SQ-29548  | BML-RA103   | 1 mg          |

## APOPTOSIS & NECROSIS

### GFP-CERTIFIED® Apoptosis/Necrosis Detection Kit (ENZ-51002)










GFP-compatible multiplex assay for distinguishing between healthy, early apoptotic, late apoptotic and necrotic cells.

- True multiplexing capabilities with GFP and other green fluorescent probes
- Optimized for both fluorescence microscopy and flow cytometry applications
- Suitable for death pathway analysis and drug/toxin studies

### NUCLEAR-ID® Blue/Red Cell Viability Reagent (GFP-CERTIFIED®) (ENZ-53005)

A mixture of a blue fluorescent cell-permeable nucleic acid dye and a red fluorescent cell-impermeable nucleic acid dye that is suited for staining dead nuclei.

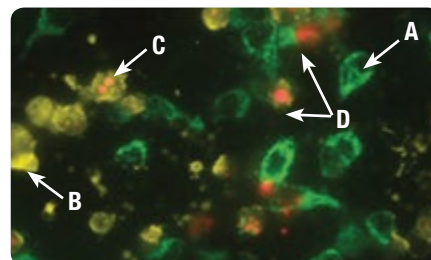
- Optimized for use in demanding imaging applications, such as confocal microscopy, flow cytometry and HCS, where consistency and reproducibility are required
- Validated for multiplexed analysis allowing for detection of inflammatory mediators and cell viability

| Assay Kits                                     |           |   |            |  |
|--|-----------|---|------------|--|
| Product Name                                   | Product # | Application   | Size       |  |
| NUCLEAR-ID® Green Chromatin Condensation Assay | ENZ-51021 |    | 200 Assays |  |
| NUCLEAR-ID® Blue/Green Cell Viability Reagent  | ENZ-53004 |     | 200 Assays |  |
| NUCLEAR-ID® Blue/Red Cell Viability Reagent    | ENZ-53005 |     | 200 Assays |  |
| NUCLEAR-ID® Red/Green Cell Viability Reagent   | ENZ-53006 |     | 200 Assays |  |



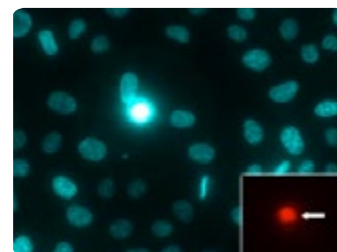
  
 Microscopy   Flow Cytometry   Microplate

### Detect Cell Death in GFP Expressing Cell Line



The Apoptosis Detection Reagent (Gold) and Necrosis Detection Reagent (Red) specifically detect cell state with clear spectral separation from mitochondria-associated GFP signal (Green).

### Convenient GFP-Compatible Viability Dye



NUCLEAR-ID® Blue/Red dye is detected as only blue-stained nuclei in live cells and fluorescent-red nuclei in dead cells (inset, arrow).

## PRO-INFLAMMATORY MARKERS

Control of pro-inflammatory molecules is critical in preventing cytotoxic effects on skin cells. Enzo provides high sensitivity ELISA kits for monitoring levels of these molecules following application of various cosmetic substances.

### IL-1 $\beta$ (human) ELISA (ADI-900-130)

- Sensitive ELISA quantifies < 1 pg/mL
- Validated for human plasma, serum and urine
- No cross-reactivity with related cytokines
- Results from up to 42 samples in duplicate in just 4 hours
- Ready-to-use reagents and pre-coated plates to reduce errors and save time

### Histamine ELISA Kit (ENZ-KIT140)

- Ultra-sensitive measurement of histamine, detecting as little as 0.03 ng/mL
- Does not require an acylation step
- Broad dynamic range suitable for a large variety of samples
- High-throughput format with results in 2 hours for up to 40 samples in duplicate

| ELISA Kits                                |              |             |
|---|--------------|-------------|
| Product Name                              | Product #    | Size        |
| Histamine ELISA Kit                       | ENZ-KIT140   | 1x96 Wells  |
| IL-1 $\beta$ (human) ELISA Kit            | ADI-900-130  | 1x96 Wells  |
| IL-1 $\beta$ (mouse) ELISA Kit            | ADI-900-132A | 1x96 Wells  |
| IL-1 $\beta$ (rat) ELISA Kit              | ADI-900-131  | 1x96 Wells  |
| IL-6 (human) ELISA Kit                    | ADI-900-033  | 1x96 Wells  |
| IL-6 (human) ELISA Kit                    | ADI-901-033  | 5x96 Wells  |
| IL-6 (mouse) ELISA Kit                    | ADI-900-045  | 1x96 Wells  |
| IL-6 (human) High Sensitivity ELISA Kit   | ENZ-KIT178   | 1x96 Wells  |
| IL-8 (human) ELISA Kit                    | ADI-900-156  | 1x96 Wells  |
| IL-8 (human) ELISA Kit                    | ADI-901-156  | 5x96 Wells  |
| IL-8 (human) ELISA Kit                    | ADI-902-156  | 10x96 Wells |
| IL-17A (human) ELISA Kit                  | ADI-900-177  | 1x96 Wells  |
| TIMP-1 ELISA Kit                          | ENZ-KIT147   | 1x96 Wells  |
| TNF- $\alpha$ (human) ELISA Kit           | ADI-900-099  | 1x96 Wells  |
| TNF- $\alpha$ (human) ELISA Kit           | ADI-901-099  | 5x96 Wells  |
| TNF- $\alpha$ (mouse) ELISA Kit           | ADI-900-047  | 1x96 Wells  |
| TNF- $\alpha$ (rat) ELISA Kit             | ADI-900-086A | 1x96 Wells  |
| TNF-R1, soluble (human) ELISA Kit         | ALX-850-047  | 1x96 Wells  |
| Antibodies                                |              |             |
| Product Name                              | Product #    | Size        |
| IL-6 (mouse), mAb (6B4 IGH 54)            | ALX-804-335  | 100 $\mu$ g |
| IL-8 receptor, Type B (human), mAb (HC 2) | ALX-804-040  | 100 $\mu$ g |
| IL-8 receptor, Type B (human), mAb (X2.7) | ALX-804-593  | 100 $\mu$ g |
| IL-17E, mAb (68C1039.2)                   | ALX-804-377  | 100 $\mu$ g |
| TNF receptor 2, pAb                       | ADI-905-593  | 1 mL        |
| TNF- $\alpha$ (human), mAb (TNF-D)        | ALX-804-199  | 100 $\mu$ g |
| TNF- $\alpha$ , mAb (1)                   | ADI-CSA-806  | 100 $\mu$ g |
| TNF- $\alpha$ , pAb                       | ALX-210-335  | 100 $\mu$ g |
| TNF-R1 (human), mAb (H398)                | ALX-804-200  | 100 $\mu$ g |
| TNF-R2, mAb (80M2)                        | ALX-804-450  | 100 $\mu$ g |

## DNA DAMAGE

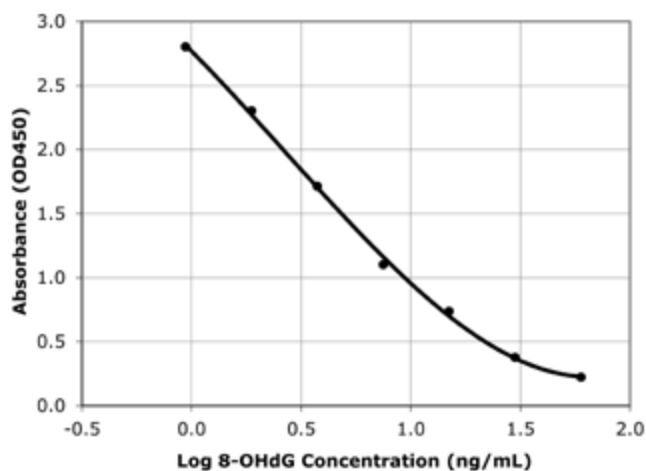
### ***Rapidly monitor DNA destruction arising from cancer, apoptosis and oxidative stress using the DNA Damage ELISA kit***

#### **DNA Damage ELISA KIT**

The DNA Damage ELISA is a fast and sensitive immunoassay providing results in less than 2.5 hours. Quantitation of 8-hydroxy-2'-deoxyguanosine (8-OHdG) in urine, serum, and saliva samples is performed in a convenient 96-well plate format using a colorimetric substrate. 8-OHdG is a frequently-used critical biomarker of oxidative stress and carcinogenesis.

- Quantify levels < 1 ng/mL
- Validated in-house in a variety of sample matrices
- Tested in a variety of biofluids (urine, serum, and saliva)
- Convenient colorimetric 96-well plate format

#### **Typical 8-OHdG Standard Curve**



The standard curve has a range of 0.94 – 60 ng/mL.

| Product Name         | Product #   | Size       |
|----------------------|-------------|------------|
| DNA Damage ELISA Kit | ADI-EKS-350 | 1x96 Wells |

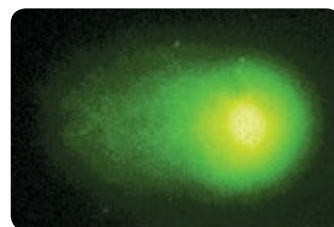
## COMET SCGE ASSAY KIT

### ***Sensitive and versatile method for measuring single- and double-strand DNA breaks in individual cells***

Exposure of cells to oxidative and environmental stresses frequently results in the breakdown or oxidation of genomic DNA. Assays to evaluate the integrity of genomic DNA, or to assess the presence of oxidized DNA are frequently used as a means of verifying the onset of apoptosis or DNA damage. The Comet SCGE Assay measures DNA damage by fluorescently detecting the integrity of DNA liberated from cells embedded in low melting point agarose. Upon electrophoresis, fragmented DNA produces a characteristic “comet” shaped tail as small DNA fragments migrate in the gel more rapidly than in-tact genomic DNA.

The Comet SCGE Assay is a fast and simple electrophoresis method to detect and quantitate DNA fragmentation in cells associated with DNA damage and apoptosis. A unique nucleic acid stain provides improved sensitivity for DNA visualization compared to ethidium bromide.

- Comet Slides allow direct application of sample without pretreatment
- Shorter assay time allows for higher throughput sample analysis
- Hydrophobic barrier allows sample treatment with DNA repair enzymes
- Unique nucleic acid stain provides improved sensitivity for DNA visualization compared to ethidium bromide

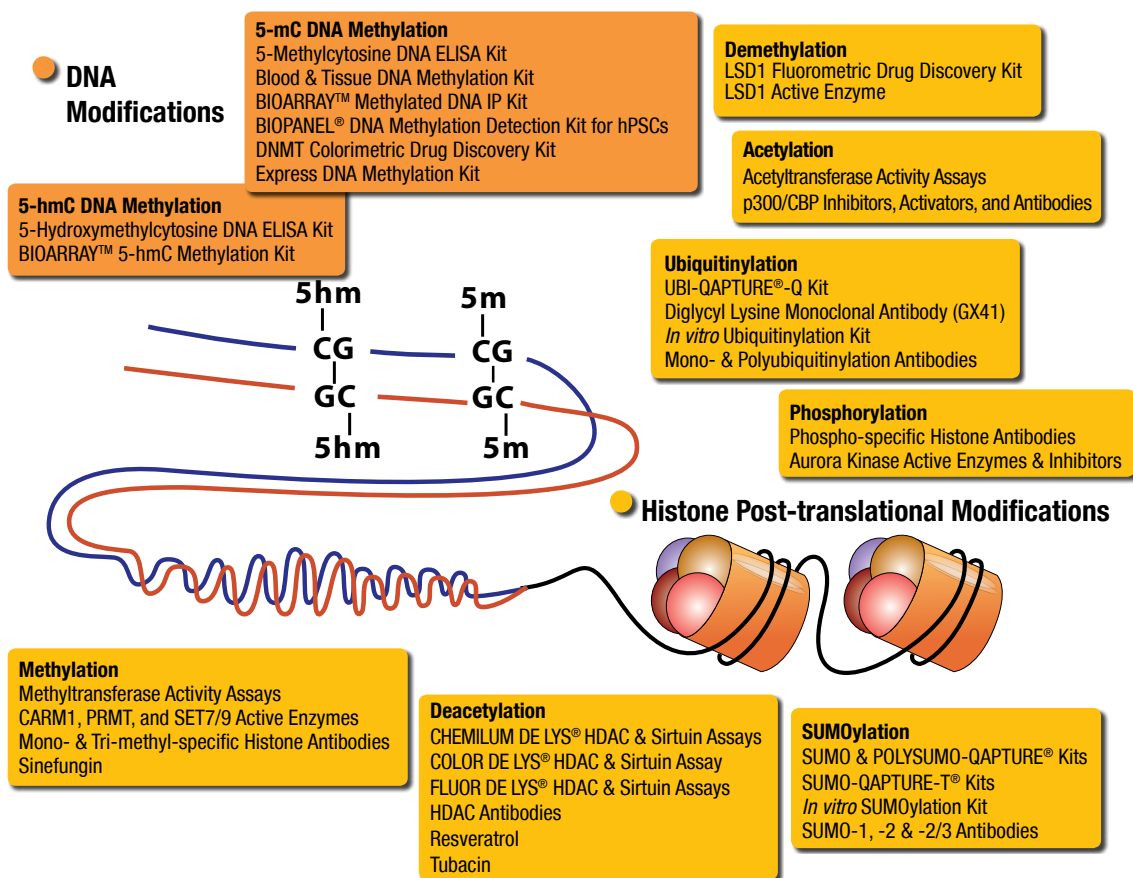


Positive control with DNA from cells treated with 100 mM H<sub>2</sub>O<sub>2</sub> for 20 minutes at 4°C.

| Product Name         | Product #   | Size     |
|----------------------|-------------|----------|
| Comet SCGE Assay Kit | ADI-900-166 | 50 Tests |

## GENE DEREGULATION

Epigenetic control of skin development and regeneration is an active area of research. Diet and exposure to environmental stresses are two significant factors in determining a change in epigenetic patterns. As such, personal care researchers are targeting epigenetic mechanisms as they formulate new products for the anti-aging market. Enzo offers a complete selection of active enzymes, antibodies and sensitivity detection assays for understanding epigenetic changes involving DNA methylation and post-translational modifications.



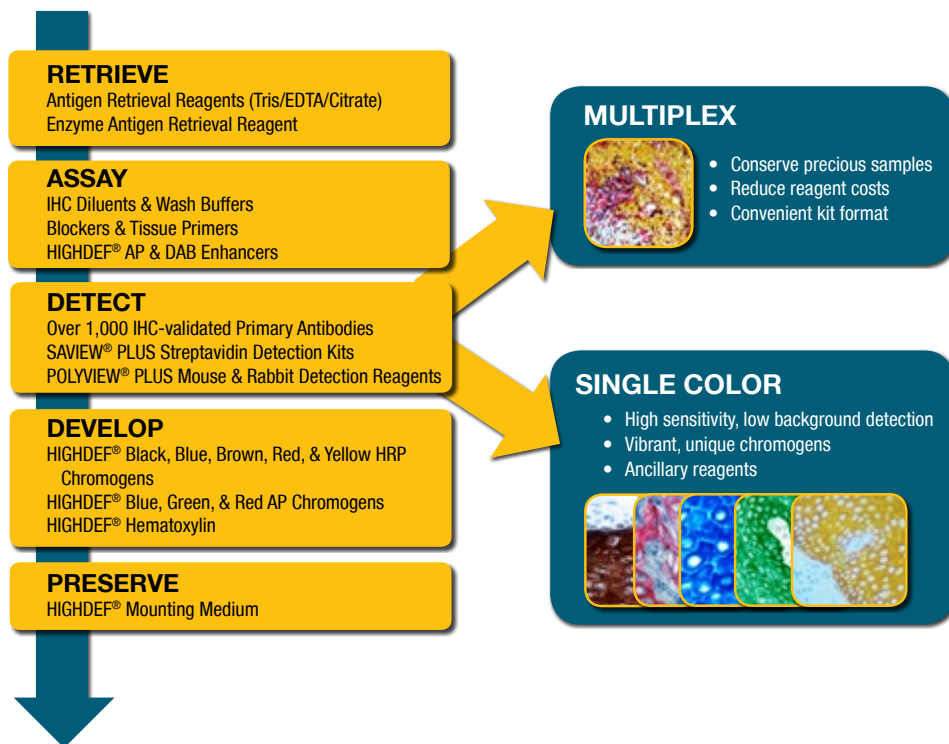
Need more information?  
Download our **EPIGENETICS** Brochures



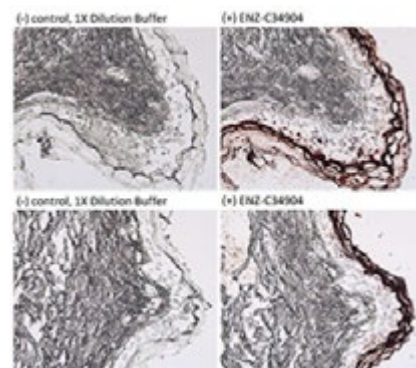
# IMMUNOHISTOCHEMISTRY FOR SKIN BIOMARKERS

## Providing the Whole Assay: Retrieval to Mounting

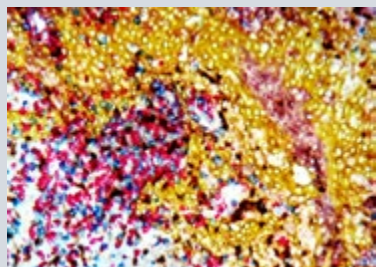
There is no need to go anywhere else for the antibodies, ancillary reagents, or other materials you need for your IHC process. Enzo Life Sciences offers a complete set of tools to help improve the quality and efficiency of your IHC process from start to finish.



### Cytokeratin 1 (human), mAb (34βB4) (ENZ-C34904)



Immunohistochemistry analysis of human skin tissue using Cytokeratin 1 (human), mAb (34βB4) at a dilution of 1:20.



Formalin-fixed paraffin-embedded human tonsil stained with pre-diluted mouse CD68 antibody labeled with HIGHDEF® IHC chromogen substrate (DAB, HC) (ADI-950-211, brown), pre-diluted rabbit lambda light-chain antibody labeled with HIGHDEF® Blue IHC chromogen (AP) (ADI-950-150, blue), pre-diluted rabbit kappa light-chain antibody labeled with HIGHDEF® Red IHC chromogen (AP, plus) (ADI-950-141, red), high molecular weight cytokeratin labeled with HIGHDEF® Yellow IHC chromogen (HRP) (ADI-950-170, yellow), and methyl green nuclear counterstain.

- Widest panel of color choices
- Unique yellow HRP chromogen
- Increased flexibility with HRP & AP options
- High definition color development

Need more information?  
Download our **IHC** Brochure

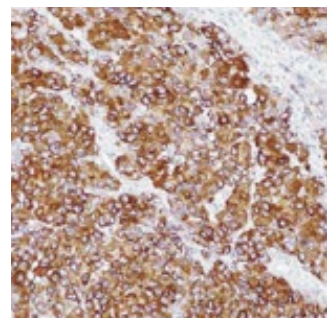




Our growing list of over 1,000 IHC-validated antibodies includes those for the detection of key skin cell markers, mediators of cell signaling and cell death, oxidative stress, heat shock proteins, proteasomes, and more. We know IHC optimization can be a time- and resource-consuming endeavor. You can back any of our antibodies with our Worry-Free Antibody Trial Program, even those not validated for IHC. Visit [www.enzolifesciences.com/abtrial](http://www.enzolifesciences.com/abtrial) for details.

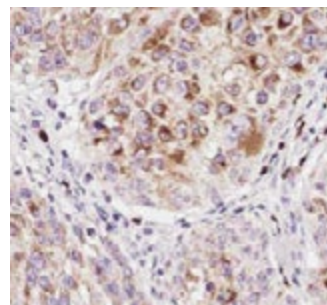
| Antibodies   |                |                     |
|--|----------------|---------------------|
| Product Name   | Product #      | Size                |
| Aggrecan, mAb (5C5)  | ALX-803-311    | 100 µL              |
| Aggrecan, mAb (5D3)  | ALX-803-313    | 100 µL              |
| CARMA1, pAb (AL220)  | ALX-804-848    | 100 µg              |
| Caspase-2, mAb (10C6)                                      | ALX-804-356    | 100 µg              |
| Caspase-3 (active), pAb                                    | BML-SA320      | 100 µL              |
| Caspase-3, mAb (31A1067)                                   | ALX-210-806    | 100 µg              |
| Caspase-3, pAb   | ADI-AAP-113    | 50 µg, 200 µg       |
| Caspase-7, mAb (10-1-62)                                   | ADI-AAM-137    | 25 µg, 100 µg       |
| Caspase-7, mAb (7-1-11)                                    | ADI-AAM-127    | 100 µg              |
| Caspase-9, mAb (2-22)                                      | ADI-AAM-139    | 100 µg              |
| Caspase-9, pAb   | ADI-AAP-149    | 25 µg, 100 µg       |
| CD44std (human), mAb (SFF-304)                             | ALX-801-089    | 100 µg              |
| CD44var(v10) (human), mAb (VFF-14)                         | ALX-801-084    | 100 µg              |
| CD44var(v3-v10) (human), pAb                               | ALX-210-234    | 1 mg                |
| Collagen type I, mAb (5D8)                                 | BPD-CSI-008-01 | 200 µg, 1 mg        |
| Collagen type III, mAb (1E7-D7/Col3)                       | BPD-CSI-007-01 | 200 µg              |
| Collagen type XII, mAb (378D5)                             | ALX-803-309    | 100 µL              |
| Cytokeratin 1 (human), mAb (34βB4) (Ready-to-Use)          | ENZ-30904      | 6 mL                |
| Cytokeratin 1 (human), mAb (34βB4)                         | ENZ-C34904     | 0.5 mL              |
| Cytokeratin 1/5/10/14 (human), mAb (34βE12)                | ENZ-C34903     | 0.5 mL              |
| Cytokeratin 1/5/10/14 (human), mAb (34βE12) (Ready-to-Use) | ENZ-30903      | 6 mL                |
| Cytokeratin 3, mAb (AE5)                                   | BML-CB10218    | 50 µg               |
| Cytokeratin 7 (human), cAb (R17-S)                         | ALX-810-218    | 1 mL                |
| Cytokeratin 8 (human), mAb (35βH11)                        | ENZ-C34902     | 0.5 mL              |
| Cytokeratin 8 (human), mAb (35βH11) (Ready-to-Use)         | ENZ-30902      | 6 mL                |
| Glutathione Peroxidase mAb (GPX-347)                       | ADI-SPA-541    | 100 µg              |
| Glutathione S-transferase-Pi mAb (USal-hGST-Pi-McAb-1)     | ALX-804-510    | 100 µg              |
| HDAC1 pAb  | ALX-804-599    | 200 µg              |
| HDAC2 pAb  | BML-SA401      | 100 µg              |
| HSP47, mAb (M16.10A1)                                      | ADI-SPA-470    | 50 µg, 200 µg, 1 mg |
| IL-1 receptor type II (human), mAb (MNC2)                  | ALX-804-461    | 100 µg              |
| IL-21R (ED), pAb   | ADI-905-308    | 100 µg              |
| IL-23 receptor, pAb  | ALX-210-636    | 100 µg              |
| IL-33, mAb (Nessy-1) (biotin conjugate)                    | ALX-804-840B   | 100 µg              |
| IL-33, mAb (Nessy-1) (preservative free)                   | ALX-804-840PF  | 100 µg              |

### MART-1 (human), cAb (A19-P)(ALX-810-217)



Formalin-fixed and paraffin-embedded human melanoma (4 µm) stained with cAb to MART-1 (human) (Prod. No. ALX-810-217) shows positive cytoplasmic immunostaining of tumor cells. Kindly performed and provided by Katarína Poliaková, MD and Lubomír Straka, MD, Ph.D. from Clinical Pathology Presov, Ltd., Presov, Slovakia.

### Melanosome (human), cAb (P14-V)(A19-P) (ALX-810-216)



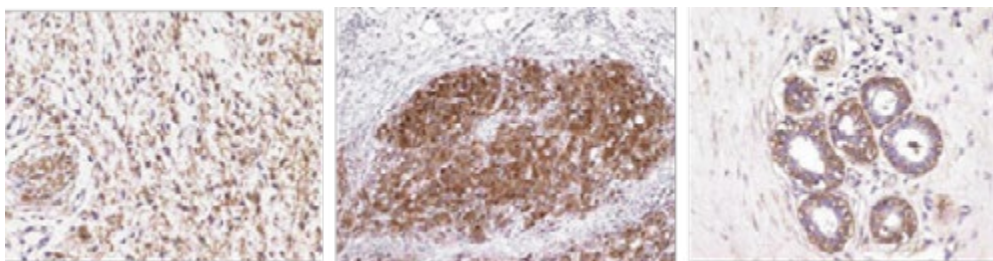
Formalin-fixed and paraffin-embedded human melanoma skin cancer tissue (4 µm) stained with cAb to melanosome (human) (Prod. No. ALX-810-216) shows specific positive immunostaining. Kindly performed and provided by Katarína Poliaková, MD and Lubomír Straka, MD, Ph.D. from Clinical Pathology Presov, Ltd., Presov, Slovakia.



| Antibodies  |               |               |
|---|---------------|---------------|
| Product Name  | Product #     | Size          |
| IL-33, mAb (Nessy-1)                                | ALX-804-840   | 100 µg        |
| IL-8 receptor, Type B (human), mAb (HC 2)           | ALX-804-040   | 50 µL         |
| LOXL4 (human), pAb                                  | ALX-215-067   | 100 µL, 1mL   |
| MART-1 (human), cAb (A19-P)                         | ALX-810-217   | 6 mL          |
| Melanoma marker (human), mAb (HMB45) (Ready-to-Use) | ENZ-30930     | 0.5 mL        |
| Melanoma marker (human), mAb (HMB45)                | ENZ-C34930    | 1 mL          |
| Melanosome (human), cAb (P14-V)                     | ALX-810-216   | 1 mL          |
| MMP-1, pAb  | ADI-905-472   | 100 µg        |
| MMP-2 (CT), pAb                                     | BML-SA616     | 500 µL        |
| MMP-9 (CT), pAb                                     | BML-SA620     | 100 µg        |
| MMP-9, pAb  | ALX-210-756   | 50 µL         |
| MMP-12 (hinge region), pAb                          | BML-SA669     | 50 µg         |
| NLRP1/NALP1 (human) (CT), pAb (Bur 242)             | ALX-210-018   | 100 µg        |
| NLRP1/NALP1 (human) (NT), pAb (Bur 241)             | ALX-804-803   | 100 µg        |
| p53 DINP1 SIP, pAb                                  | ADI-905-300   | 50 µg, 200 µg |
| p53, mAb (PAb122)                                   | ADI-KAM-CC002 | 250 µg        |
| p53, pAb  | ADI-KAP-CC030 | 1 mL          |
| p53, pAb  | ADI-905-510   | 100 µg        |
| RAIDD, pAb  | ADI-905-244   | 500 µL        |
| S100 (human), cAb (D28-E)                           | ALX-810-220   | 100 µL        |
| SIRT1 (human), pAb                                  | BML-SA427     | 100 µL        |
| SIRT3 pAb   | BML-SA463     | 100 µg        |
| Smac/DIABLO (CT), pAb                               | ADI-905-684   | 100 µg        |
| Smac/DIABLO (human), pAb                            | ALX-804-366   | 100 µL        |
| Smac/DIABLO, pAb                                    | BML-SA101     | 100 µL        |

### S100 (human), cAb (D28-E)(ALX-810-220)

S100 is useful in understanding the origins of various skin tumors and in assessing the diagnosis and prognosis of a variety of skin lesions.



Formalin-fixed and paraffin-embedded human neurofibroma tissue (left; 4 µm), skin melanoma tissue (middle; 4 µm), and human breast tissue (right; 4 µm) stained with cAb to S100 (human) (Prod. No. ALX-810-220) show: Left - positive immunostaining of tumor cells and peripheral nerve; Middle - strong positive immunostaining of melanoma cells; Right - strong positive immunostaining of the outer myoepithelial cell component in mammary lobule. Kindly performed and provided by Katarína Poliaková, MD and Lubomír Straka, MD, Ph.D. from Clinical Pathology Presov, Ltd., Presov, Slovakia.

# ENDOCRINE DISRUPTION

## HORMONES

### Rapid Identification of Endocrine Disruptors Using High-Sensitivity ELISAs

Studies indicate that environmental chemicals, such as those found in personal care products and packaging, can interfere with the production of hormones in humans and other animals. Enzo Life Sciences provides high-sensitivity ELISA kits and complementary reagents for researchers looking to assess the potential endocrine disruptive effects of substances.

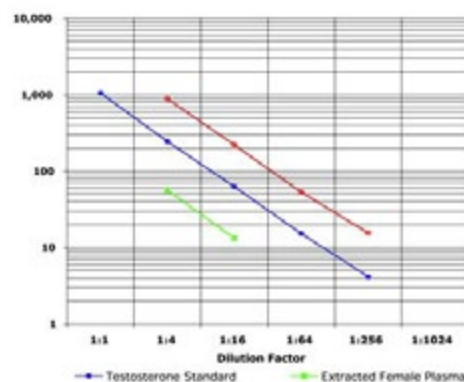
#### Testosterone High Sensitivity ELISA Kit (ADI-900-176)

A highly sensitive competitive immunoassay for the measurement of testosterone in human serum and plasma samples.

- Highest sensitivity ELISA on the market (3.9 – 1000 pg/mL)
- Validated for use with human samples types (plasma, serum, and urine)
- Includes stable color-coded reagents to reduce user errors

| Hormone ELISA Kits                      |             |            |
|---|-------------|------------|
| Product Name                            | Product #   | Size       |
| Testosterone ELISA Kit                  | ADI-900-065 | 1x96 Wells |
|   | ADI-901-065 | 5x96 Wells |
| Testosterone High Sensitivity ELISA Kit | ADI-900-176 | 1x96 Wells |
|   | ADI-901-176 | 5x96 Wells |

#### High Specificity Detection of Testosterone



Dose-response curves from extracted male serum and extracted female plasma diluted into assay buffer were compared to the Testosterone standard curve. The parallel response indicates that the standard effectively mimics the native protein.

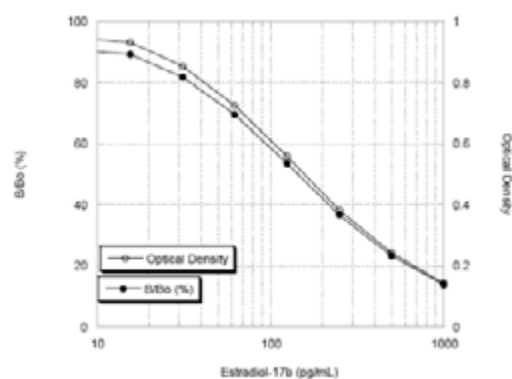
#### 17β-Estradiol High Sensitivity ELISA Kit (ADI-900-174)

The 17β-Estradiol (serum/plasma) ELISA is a colorimetric competitive enzyme immunoassay kit for rapid detection of this analyte in human samples.

- High sensitivity — detects down to 14.0 pg/mL
- Validated for use with human samples types (plasma and serum)
- Fast, reliable results from up to 37 samples in duplicate in just 3 hours

| Hormone ELISA Kits                       |             |            |
|--|-------------|------------|
| Product Name                             | Product #   | Size       |
| 17β-Estradiol ELISA Kit                  | ADI-900-008 | 1x96 Wells |
|  | ADI-901-008 | 5x96 Wells |
| 17β-Estradiol High Sensitivity ELISA Kit | ADI-900-174 | 1x96 Wells |
|  | ADI-901-174 | 5x96 Wells |

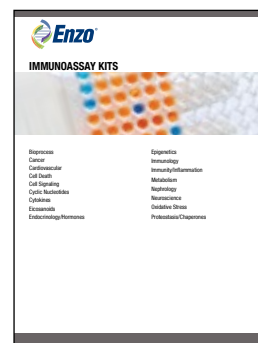
#### Detect Picogram levels of 17β-Estradiol



Plot of the Percent Bound (B/B<sub>0</sub>) versus concentration of 17β-Estradiol for the standards. The amount of signal is indirectly proportional to the amount of 17β-Estradiol in the sample.

| Hormone ELISA Kits   |                              |                          |
|--|------------------------------|--------------------------|
| Product Name   | Product #                    | Size                     |
| 25(OH) Vitamin D ELISA Kit   | ADI-900-215                  | 1x96 Wells               |
| Arg <sup>8</sup> -Vasopressin ELISA Kit                                  | ADI-900-017A<br>ADI-901-017A | 1x96 Wells<br>5x96 Wells |
| Big Endothelin-1 (human), ELISA Kit<br>Big Endothelin-1 (rat), ELISA Kit | ADI-900-022<br>ADI-900-073   | 1x96 Wells<br>1x96 Wells |
| Corticosterone ELISA Kit   | ADI-900-097<br>ADI-901-097   | 1x96 Wells<br>5x96 Wells |
| Cortisol ELISA Kit   | ADI-900-071<br>ADI-901-071   | 1x96 Wells<br>5x96 Wells |
| DHEA ELISA Kit   | ADI-900-093<br>ADI-901-093   | 1x96 Wells<br>5x96 Wells |
| Endothelin-1 ELISA Kit   | ADI-900-020A                 | 1x96 Wells               |
| Estriol ELISA Kit  | ADI-900-100                  | 1x96 Wells               |
| Gastrin I (human), ELISA Kit<br>Gastrin I (rat), ELISA Kit               | ADI-900-026<br>ADI-900-149   | 1x96 Wells<br>1x96 Wells |
| Leptin (human), ELISA Kit<br>Leptin (mouse), ELISA Kit                   | ADI-900-028A<br>ADI-900-019A | 1x96 Wells<br>1x96 Wells |
| Leptin (rat), ELISA Kit  | ADI-900-015A                 | 1x96 Wells               |
| Oxytocin ELISA Kit   | ADI-900-153A<br>ADI-901-153A | 1x96 Wells<br>5x96 Wells |
| Progesterone ELISA Kit   | ADI-900-011<br>ADI-901-011   | 1x96 Wells<br>5x96 Wells |
| Serotonin ELISA Kit  | ADI-900-175                  | 1x96 Wells               |
| Substance P ELISA Kit  | ADI-900-018<br>ADI-901-018   | 1x96 Wells<br>5x96 Wells |

Need more information?  
Download our **IMMUNOASSAY KITS**  
Product Guide



# ENDOCRINE DISRUPTION

| Hormone Antibodies   |                  |        |
|--|------------------|--------|
| Product Name   | Product #        | Size   |
| [pTyr <sup>1150</sup> /Tyr <sup>1151</sup> ]Insulin receptor, mAb (10C3)                                   | ADI-905-645      | 100 µg |
| [pTyr <sup>1158/1162/1163</sup> ]Insulin receptor and [pTyr <sup>1131/1135/1136</sup> ]IGF-1 Receptor, pAb | BML-SA392        | 20 µL  |
| [pTyr <sup>1322</sup> ]Insulin receptor, mAb (21G12)   | ADI-905-646      | 100 µg |
| Angiotensin II receptor AT1, pAb   | ADI-905-743      | 100 µg |
| Angiotensin II receptor AT1, pAb (DY-682 conjugate)  | ADI-905-813      | 100 µg |
| Angiotensin II receptor AT1, pAb (DY-800 conjugate)  | ADI-905-814      | 100 µg |
| Angiotensin II receptor AT2, pAb   | ADI-905-746      | 100 µg |
| Angiotensin II receptor AT2, pAb (DY-682 conjugate)  | ADI-905-815      | 100 µg |
| AT1 receptor (extracellular), pAb  | BML-SA608        | 50 µL  |
| AT2 receptor (extracellular), pAb  | BML-SA609        | 50 µL  |
| Atrial natriuretic peptide, mAb (23/1)   | BML-AA6244       | 100 µL |
| Atrial natriuretic peptide, pAb  | BML-AA1105       | 25 µL  |
| Brain natriuretic peptide, pAb   | BML-BA1117       | 25 µL  |
| Calcitonin gene related peptide, mAb (CD8)   | BML-CA1135       | 25 µL  |
| Calcitonin gene related peptide, pAb   | BML-CA1134       | 25 µL  |
| Calcitonin gene related peptide, pAb   | BML-CA1137       | 25 µL  |
| Endothelin A receptor, pAb   | ALX-210-507A     | 250 µg |
| Endothelin B receptor, pAb   | ALX-210-506A     | 250 µg |
| Endothelin receptor A (rat), pAb   | BML-SA576        | 50 µL  |
| Endothelin receptor B (rat), pAb   | BML-SA577        | 50 µL  |
| Endothelin receptor ETA, pAb   | ADI-905-790      | 100 µg |
| Endothelin receptor ETB, pAb   | ADI-905-791      | 100 µg |
| Endothelin-1, mAb (TR.ET.48.5)   | ALX-803-001      | 100 µL |
| Estrogen receptor $\alpha$ , mAb (33)  | ALX-803-004      | 50 µg  |
| Estrogen receptor $\alpha$ , mAb (C-542)   | ADI-SRA-1010     | 50 µg  |
| Estrogen receptor $\alpha$ , pAb   | ALX-210-201      | 50 µg  |
| Estrogen receptor $\beta$ , pAb  | ALX-210-132      | 100 µg |
| Estrogen receptor $\beta$ , pAb  | ALX-210-135      | 50 µg  |
| Estrogen receptor $\beta$ , pAb  | ALX-210-178      | 50 µg  |
| Estrogen receptor $\beta$ , pAb  | ALX-210-180      | 50 µg  |
| Estrogen receptor, mAb (h-151)   | ADI-SRA-1000     | 50 µg  |
| Galanin, pAb   | BML-GA1161       | 25 µL  |
| Gastrin releasing peptide, pAb   | BML-GA1166       | 25 µL  |
| Gastrin, pAb   | BML-GA1167       | 25 µL  |
| Ghrelin (human), mAb (45)  | BPD-ABS-050-45   | 200 µg |
| Ghrelin receptor GHS-R, pAb  | ADI-905-792      | 100 µg |
| Ghrelin, mAb (121)   | BPD-ABS-052-121  | 200 µg |
| Ghrelin, mAb (121) (biotin conjugate)  | BPD-ABS-052-121B | 50 µg  |
| Glucagon like peptide 1, pAb   | BML-GA1176       | 25 µL  |
| Glucagon like peptide 2, pAb   | BML-GA1179       | 25 µL  |

| Hormone Antibodies                                |                 |        |
|---|-----------------|--------|
| Product Name                                      | Product #       | Size   |
| Glucagon, mAb (08)                                | BPD-ABS-058-08  | 200 µg |
| Glucagon, mAb (08) (biotin conjugate)             | BPD-ABS-058-08B | 50 µg  |
| Glucagon, mAb (09)                                | BPD-ABS-061-09  | 200 µg |
| Glucagon, pAb                                     | BML-GA1181      | 25 µL  |
| Glucocorticoid receptor α, pAb                    | ALX-210-277     | 50 µg  |
| Glucocorticoid receptor β (human), pAb            | ALX-210-154     | 50 µL  |
| Glucocorticoid receptor, mAb (BugR2)              | ALX-803-006     | 100 µg |
| Glucose-dependent insulinotropic peptide, pAb     | BML-GA1173      | 25 µL  |
| Gonadotropin releasing hormone, mAb (BML)         | BML-GZ1092      | 25 µL  |
| Gonadotropin releasing hormone, pAb               | BML-GZ1110      | 25 µL  |
| Insulin receptor β subunit, mAb (CT-3)            | BML-SA432       | 100 µg |
| Insulin receptor β, mAb (C18C4)                   | ADI-905-683     | 100 µg |
| Insulin receptor, mAb (9H4)                       | ADI-905-647     | 100 µg |
| Leptin receptor, pAb                              | ALX-210-168     | 100 µL |
| Leptin, pAb                                       | ALX-210-105     | 50 µL  |
| Leptin, pAb                                       | BML-SA268       | 100 µg |
| Nesfatin-1, mAb (Nesfaty-1)                       | ALX-804-854     | 100 µg |
| Nesfatin-1, pAb (AT120)                           | ALX-210-944     | 100 µg |
| Orexin receptor 1, pAb                            | BML-SA646       | 50 µL  |
| Orexin receptor 2, pAb                            | BML-SA647       | 50 µL  |
| Progesterone receptor, mAb (aPR6)                 | ALX-803-010     | 100 µg |
| Progesterone, mAb (111.2C7.3)                     | ADI-905-010     | 100 µL |
| Prolactin receptor (rat), mAb (T6)                | ALX-803-003     | 100 µg |
| Prolactin receptor, mAb (U5)                      | ALX-803-005     | 100 µg |
| Serotonin receptor 5HT1A pAb                      | ADI-905-741     | 100 µg |
| Serotonin, pAb                                    | BML-SZ1011      | 25 µL  |
| Serotonin, pAb                                    | BML-SZ1021      | 100 µL |
| Somatostatin 28, pAb                              | BML-SA1268      | 100 µL |
| Somatostatin receptor type 3, pAb                 | BML-SA593       | 50 µL  |
| Somatostatin receptor type 4 (rat), pAb           | BML-SA594       | 200 µL |
| Somatostatin, pAb                                 | BML-SZ1114      | 100 µL |
| Somatostatin, pAb                                 | BML-SZ1116      | 25 µL  |
| Substance P, pAb                                  | BML-SA1270      | 25 µL  |
| Substance P, pAb                                  | BML-SZ1062      | 100 µL |
| Thyroid Peroxidase (human), mAb (TPO 47)          | ALX-BC-6010-S   | 1 mL   |
| Viable rat beta cells and RIN cells, mAb (K14D10) | ALX-803-052     | 100 µg |
| Vitamin D3 receptor, mAb (9A7)                    | ALX-804-021     | 100 µg |
| WIPI-1 (human), pAb (AT129)                       | ALX-210-954     | 50 µg  |
| WIPI-1, pAb (AT130)                               | ALX-210-955     | 100 µg |

## BIOCHEMICALS

| Biochemicals                                |             |                     |
|---|-------------|---------------------|
| Product Name                                | Product #   | Size                |
| (R,S)-Equol                                 | ALX-385-032 | 5 mg, 25 mg         |
| (Z)-4-Hydroxytamoxifen                      | ALX-550-361 | 1 mg, 5 mg          |
| 1 $\alpha$ ,25-Dihydroxyvitamin D3          | BML-DM200   | 50 $\mu$ g, 1 mg    |
| 2-[1-(4-Piperonyl)piperazinyl]benzothiazole | ALX-550-171 | 5 mg                |
| 20-Hydroxyecdysone                          | ALX-370-012 | 5 mg, 10 mg, 50 mg  |
| 24(R),25-Dihydroxyvitamin D3                | BML-DM300   | 50 $\mu$ g          |
| 25-Hydroxyvitamin D2                        | BML-DM101   | 1 mg                |
| 25-Hydroxyvitamin D3                        | BML-DM100   | 1 mg, 5 mg          |
| 2-Methoxyestradiol                          | BML-S540    | 10 mg, 50 mg        |
| 6 $\alpha$ -Fluorotestosterone              | BML-S250    | 5 mg, 25 mg         |
| 8-Isopentenylnaringenin                     | ALX-385-025 | 5 mg                |
| Alfacalcidol                                | ALX-460-036 | 1 mg                |
| Bicalutamide                                | ALX-270-476 | 100 mg, 500 mg, 1 g |
| Compound A                                  | ALX-550-516 | 5 mg                |
| Corticosterone                              | ALX-370-007 | 50 mg               |
| Coumestrol                                  | BML-S180    | 5 mg, 25 mg         |
| Dihydroergocristine . mesylate              | BML-NS108   | 200 mg, 1 g         |
| DPN   | BML-GR248   | 10 mg, 50 mg        |
| Ferutinin (high purity)                     | ALX-350-098 | 1 mg, 5 mg, 10 mg   |
| Finasteride                                 | ALX-270-491 | 100 mg, 500 mg      |
| Formononetin (high purity)                  | ALX-270-312 | 5 mg                |
| HSD Inhibitor 23                            | ALX-270-474 | 1 mg                |
| Isoxanthohumol                              | ALX-350-279 | 1 mg                |
| Makisterone A                               | ALX-370-013 | 1 mg                |
| Megestrol acetate                           | BML-S515    | 1 g                 |
| Melengestrol acetate                        | BML-S520    | 200 mg, 1 g         |
| Mifepristone                                | BML-S510    | 25 mg, 100 mg       |
| ORG-12962                                   | BML-NS705   | 10 mg, 50 mg        |
| Ponasterone A                               | ALX-370-014 | 5 mg                |
| Raloxifene . HCl                            | BML-GR243   | 50 mg, 500 mg       |
| Tamoxifen . citrate                         | ALX-550-095 | 1 g                 |
| Vitamin D2                                  | ALX-460-025 | 1 g                 |
| Vitamin D3                                  | ALX-460-026 | 1 g, 5 g            |
| WAY-200070                                  | BML-GR247   | 10 mg, 50 mg        |

# PIGMENTATION

## Sensitive Detection for Modulators of Pigmentation

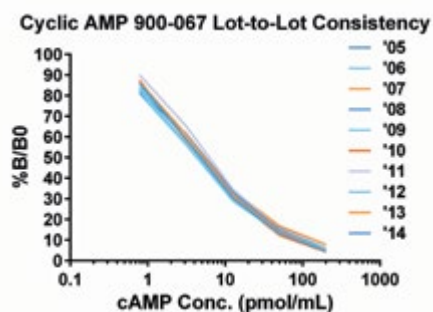
Binding of Melanocyte-Stimulating Hormone (MSH) to the melanocortin 1 receptor (MC1R, a G protein-coupled receptor) on melanocytes triggers pro-differentiation signals mediated by the second messenger cyclic AMP (cAMP). cAMP-mediated signals inhibit UV-induced apoptosis and promote melanin synthesis via transcriptional activation of pigment-regulating enzymes such as tyrosinase. Modulation of cAMP levels by pharmacological induction of adenylate cyclases, or inhibition of phosphodiesterases, hold promise as UV-independent mechanisms of increasing natural production of melanin and protection from UV damage to skin.

### Cyclic Nucleotide Analysis

Enzo Life Sciences offers the most sensitive and complete colorimetric ELISA kits for quantification of intracellular and/or extracellular cAMP or cGMP in a variety of sample types.

- Sensitive ELISA with optional acetylation protocol increases sensitivity 10-fold (0.039 pmol/mL)
- Regularly cited in peer-reviewed publications
- Simple, efficient and well-established sample handling protocols
- Reliable manufacturing yields consistent lot-to-lot performance

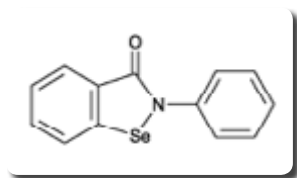
## Measure cAMP, a Regulator of Melanin Synthesis



Graph demonstrates the robust and reproducible nature of the competitive cAMP ELISA kit (Prod. #ADI-900-067) showing standard curves from 10 lots manufactured over 6 years.

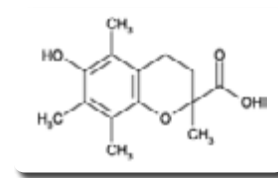
| Related Products                               |              |              |
|--|--------------|--------------|
| Product Name                                   | Product #    | Size         |
| cAMP Complete ELISA Kit                        | ADI-900-163  | 1x96 Wells   |
| cAMP Direct ELISA Kit                          | ADI-900-066  | 1x96 Wells   |
| cAMP ELISA Kit                                 | ADI-900-067  | 1x96 Wells   |
| Catalase Fluorometric Detection Kit            | ADI-907-027  | 500 Tests    |
| PKA Kinase Activity Kit                        | ADI-EKS-390A | 1x96 Wells   |
| RO-20-1724                                     | BML-EI117    | 100 mg, 1 g  |
| Rolipram                                       | BML-PD175    | 10 mg, 50 mg |
| R-(-)-Rolipram                                 | BML-PD177    | 5 mg         |
| SQ 20009                                       | BML-PD130    | 5 mg, 25 mg  |
| S-(+)-Rolipram                                 | BML-PD178    | 5 mg         |
| Zardaverine                                    | BML-PD195    | 5 mg, 25 mg  |
| PDE4A (catalytic domain) (human), recombinant  | BML-SE521    | 20 µg        |
| PDE4B2 (catalytic domain) (human), recombinant | BML-SE522    | 20 µg        |
| PDE4D (catalytic domain) (human), recombinant  | BML-SE523    | 20 µg        |

### Ebselen (ALX-270-097) - 5 mg, 25 mg



Recent findings suggest that ebselen functions as a depigmenting compound by inhibiting melanin synthesis and melanosome transfer to keratinocytes. Purity:  $\geq 98\%$  (NMR)

### Trolox® (ALX-270-267) - 100 mg



This cell-permeable vitamin E derivative decreases post-inflammatory hyperpigmentation after injury through anti-inflammatory action. Purity:  $\geq 97\%$



# SAFETY AND TOXICITY TESTING

## CYTOTOXICITY

### Whole Cell Analysis of Cytotoxic Compounds

Toxicity assessment determines whether a drug or compound induces inflammatory or cytotoxic processes such as apoptosis, cell proliferation, and oxidative stress. Enzo offers a diverse collection of assays for analysis of cell viability and cytotoxicity.

#### MITO-ID® Membrane Potential Cytotoxicity Assay Kit (ENZ-51019)

A real-time mitochondrial membrane potential assay with superior sensitivity

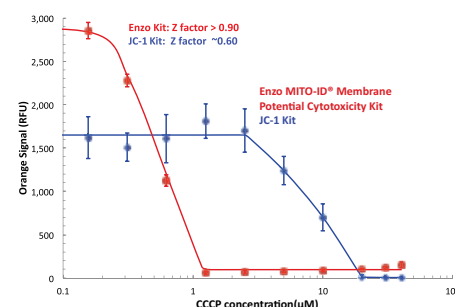
- 10x more sensitive than JC-1 with superior aqueous solubility
- Photostable dual-emission dye
- No-wash/no-medium removal
- Separate MITO-ID® Red/Green assays available for detection of mitochondrial mass
- Suitable for high-throughput applications

#### LYSO-ID® Red Cytotoxicity Kit (ENZ-51015)

A rapid, quantitative live cell cytotoxicity assay compatible with high-throughput screening (HTS)

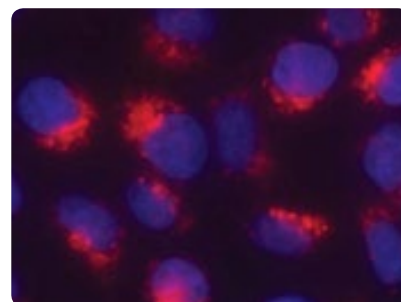
- LYSO-ID® dye rapidly partitions into cells and labels acidic organelles
- Only commercial assay available that allows for long-term cell monitoring of cytotoxic effects
- Multi-well, HTS-compatible assay with rapid 10-15 minute dye incubation

### High-throughput Analysis of Mitotoxicity














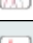
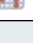








Detect mitochondrial perturbations with 10 times more sensitivity than JC-1.

### Monitor Dysfunction of Lysosomal Degradation



Drug-induced lysosome accumulation in U-2 OS cells was evaluated using Verapamil (200 μM) and LYSO-ID® Red dye.

| Assay Kits                                      |            |   |
|---|------------|---|
| Product Name                                    | Product #  | Application   |
| CYTO-ID® Autophagy Detection Kit                | ENZ-51031  |    |
| GFP-CERTIFIED® Apoptosis/Necrosis Detection Kit | ENZ-51002  |     |
| LDH Cytotoxicity WST Assay                      | ENZ-KIT157 |     |
| LYSO-ID® Red Cytotoxicity Assay                 | ENZ-51015  |     |
| MITO-ID® Membrane Potential Cytotoxicity Assay  | ENZ-51019  |    |
| MITO-ID® Membrane Potential Detection Kit       | ENZ-51018  |     |
| NUCLEAR-ID® Green Chromatin Condensation Assay  | ENZ-51021  |    |
| NUCLEAR-ID® Blue/Green Cell Viability Reagent   | ENZ-53004  |     |
| NUCLEAR-ID® Blue/Red Cell Viability Reagent     | ENZ-53005  |     |
| NUCLEAR-ID® Red/Green Cell Viability Reagent    | ENZ-53006  |     |

# NATURAL PRODUCTS AND COMPOUND SCREENING

Natural products are an unsurpassed source of chemical diversity for pharmaceutical, cosmetics, nutraceutical screening, and research. Historically, these compounds have been the most successful source of new drugs.

## SCREEN-WELL® Natural Product Library (BML-2865)

A collection of over 500 compounds of known structure in a ready-to-screen format, featuring:

- More than 500 compounds supplied in DMSO at 2.0 mg/mL
- Available in 100 µL and 500 µL sizes
- Individual compounds or custom subsets also available

| Compound Libraries            |           |                |
|-------------------------------|-----------|----------------|
| Product Name                  | Product # | Size           |
| Autophagy Library             | BML-2837  | 100 µL, 500 µL |
| Bioactive Lipid Library       | BML-2800  | 100 µL, 500 µL |
| Epigenetics Library           | BML-2836  | 100 µL, 500 µL |
| Fatty Acid Library            | BML-2803  | 100 µL, 500 µL |
| FDA Approved Library          | BML-2843  | 100 µL         |
| ICCB Known Bioactives Library | BML-2840  | 100 µL         |
| Ion Channel Library           | BML-2805  | 500 µL         |
| Natural Products Library      | BML-2865  | 100 µL, 500 µL |
| Protease Inhibitor Library    | BML-2833  | 100 µL, 500 µL |
| REDOX Library                 | BML-2835  | 100 µL, 500 µL |

| Polyphenols & Related Compounds    |             |                                  |
|------------------------------------|-------------|----------------------------------|
| Product Name                       | Product #   | Size                             |
| Delphinidin chloride (high purity) | ALX-385-028 | 10 mg                            |
| EGCG                               | ALX-270-263 | 10 mg, 50 mg                     |
| Genistein (synthetic)              | ALX-350-006 | 10 mg, 25 mg, 50 mg, 100 mg, 1 g |
| Hydroxytyrosol                     | ALX-350-404 | 10 mg, 50 mg, 250 mg             |
| Myricetin                          | ALX-385-012 | 10 mg, 50 mg                     |
| Quercetin                          | ALX-385-001 | 5 g, 25 g                        |
| Resveratrol                        | BML-FR104   | 100 mg, 500 mg                   |

| Antioxidants                                     |             |              |
|--|-------------|--------------|
| Product Name                                     | Product #   | Size         |
| Retinoic Acid, all trans                         | BML-GR100   | 500 mg, 5 g  |
| Rosmarinic Acid                                  | ALX-270-253 | 10 mg, 50 mg |
| Ubiquinone-10/Coenzyme Q                         | ALX-270-295 | 2 mg         |
| Vitamin E/DL- $\alpha$ -Tocopherol (high purity) | ALX-460-018 | 1 g          |

### ***Need a Bulk Compound?***

We offer over 2,500 biologically characterized compounds that are available in bulk upon request.

**Inquire at: [www.enzolifesciences.com/bulk](http://www.enzolifesciences.com/bulk)**



**Global Headquarters**  
**ENZO LIFE SCIENCES, INC.**  
10 Executive Blvd.  
Farmingdale, NY 11735  
Ph: 800.942.0430  
Fax: 631.694.7501  
info-usa@enzolifesciences.com

**European Sales Office**  
**ENZO LIFE SCIENCES (ELS) AG**  
Industriestrasse 17  
CH-4415 Lausen, Switzerland  
Ph: +41 61 926 8989  
Fax: +41 61 926 8979  
info-eu@enzolifesciences.com

#### **LOCAL EUROPEAN OFFICES**

**Belgium, The Netherlands & Luxembourg**  
Enzo Life Sciences BVBA  
Avenue Louise 65/Box 11  
1050 Bruxelles  
Belgium  
Ph: +32 3 466 0420  
Fax: +32 3 808 7033  
info-be@enzolifesciences.com

**France**  
Enzo Life Sciences (ELS) AG  
Branch Office Lyon  
13, avenue Albert Einstein,  
F-69100 Villeurbanne, France  
Ph: +33 472 440 655  
Fax: +33 481 680 254  
info-fr@enzolifesciences.com

**Germany**  
Enzo Life Sciences GmbH  
Basler Strasse 57a  
DE-79540 Lörrach  
Germany  
Ph: +49 7621 5500 526  
Fax: +49 7621 5500 527  
info-de@enzolifesciences.com

**UK & Ireland**  
Enzo Life Sciences (UK) Ltd.  
1 Colleton Crescent  
Exeter EX2 4DG  
Ph: 0845 601 1488 (UK customers)  
Ph: +44 1392 825900  
Fax: +44 1392 825910  
info-uk@enzolifesciences.com

For local distributors and detailed product information visit us online:  
**[www.enzolifesciences.com](http://www.enzolifesciences.com)**