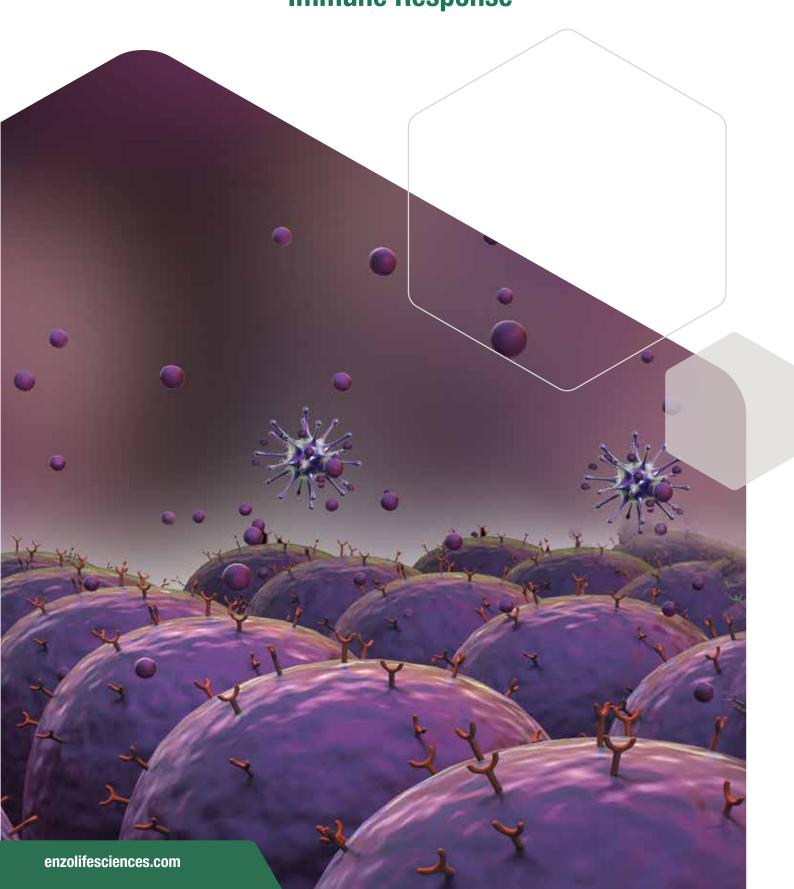
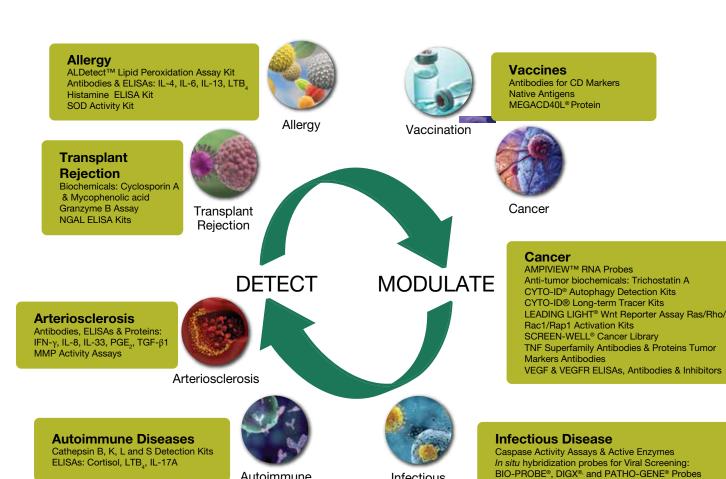


Immunology Innovative Products to Study the Immune Response



COMPREHENSIVE TOOLS FOR INTERPRETING IMMUNOLOGICAL RESPONSES

Maintaining the proper balance between immune activation and suppression is critical for the physiological function of the innate and adaptive immune system. Disruption of this equilibrium leads to immunological disorders including autoimmune diseases, hypersensitivities, and immune deficiency. Enzo provides a comprehensive portfolio of products to detect and modulate the complex processes of immunological disease.



Autoimmune

Disease

Cell Analysis & Detection

Our unique fluorescent probe-based assays and widely cited antibodies enable multiplex detection of cellular markers and key parameters of cell viability, cell signaling, death pathways, and toxicity. Our expertise in fluorescent probe chemistry and cellular analysis combined provide high-quality, flexible detection platforms for immune cell profiling and functional analysis.

Immunoassays & Antibodies

Infectious

Diseases

As a trusted manufacturer of thousands of widely cited and thoroughly validated ELISA kits and antibodies, we understand quality means delivering sensitivity, specificity, and consistency. We have over 250 immunoassays and detection kits, 3,000 antibodies, and a complete set of IHC/ ISH reagents for the analysis of immune

Small Molecule Chemistry

Toll-like Receptor Ligands: Flagellin, LPS, Lipid A, MALP-2

SCREEN-WELL® FDA Approved Drug Library

AMPIVIEW™ RNA Probes

We offer more than 3,000 biologically characterized small molecules and the industry's most diverse collection of compound libraries for predictive toxicology. Our catalog includes hundreds of redox compounds, natural products, and bioactive lipids available as ready-to-screen SCREEN-WELL® libraries or as individual and bulk compounds.

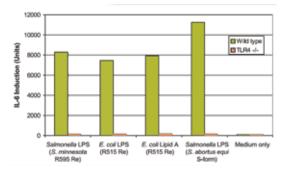
HIGH PURITY LIGANDS

Toll-like Receptors

Providing the Purest Ligands for Recognition by Pattern Recognition Receptors

| | Key Toll-like Receptor Ligands | | | | | |
|--|--------------------------------|---|-------------|--|--|--|
| | TLR | Ligands | Product # | | | |
| LPS & Lipid A Reagents | TLR 1/2 | Pam ₃ Cys-Ser-(Lys) ₄ . trihydrochloride | ALX-165-066 | | | |
| • Ultrapura 6 00 0W) | TLR 2/6 | MALP-2 | ALX-162-027 | | | |
| • <i>Ultrapure</i> (≥99.9%) | TLR 3 | Polyinosinic-polycytidylic acid. potassium salt (TLRGRADE®) (synthetic) | ALX-746-021 | | | |
| Sterile, ready-to-use liquid formulation | TLR 4 | Paclitaxel | BML-T104 | | | |
| • No hazardous handling TLR 5 TLR 7/8 | Flagellin (high purity) | ALX-522-058 | | | | |
| | R-848 | ALX-420-038 | | | | |
| Widely cited in peer-reviewed literature | TLDO | ODN 2006 (TLRGRADE®) (synthetic) | ALX-746-006 | | | |
| merature | TLR 9 | ODN 1826 (TLRGRADE®) (synthetic) | ALX-746-002 | | | |
| | TLR 11 | Profilin (Toxoplasma gondii), (recombinant) (GST-tag) | ALX-522-093 | | | |

High-Purity TLR4-activating LPS & Lipid A Preparations Do Not Activate TLR2 or other TLRs



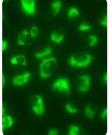
Activation of macrophages from TLR4 wild type (green) compared to TLR4 deficient mice (orange) by TLRGRADE® LPS and TLRGRADE® Lipid A.

| RELATED PRODUCTS | |
|--|-------------|
| Product Name | Product # |
| LPS from $\it E. coli$, Serotype EH100 (Ra) (TLRGRADE®) (Ready-to-Use) | ALX-581-010 |
| LPS from <i>E. coli</i> , Serotype R515 (Re) (TLRGRADE®) (Ready-to-Use) | ALX-581-007 |
| LPS from <i>E. coli</i> , Serotype 0111:B4 (TLRGRADE®) (Ready-to-Use) | ALX-581-012 |
| LPS from <i>E. coli</i> , Serotype 055:B5 (TLRGRADE®) (Ready-to-Use) | ALX-581-013 |
| LPS from S. abortus equi S-form (TLRGRADE®) (Ready-to-Use) | ALX-581-009 |
| LPS from <i>S. abortus</i> equi S-form (TLRGRADE®) (Ready-to-Use) (Biotin) | ALX-581-150 |
| LPS from $\textit{S. minnesota}$ R595 (Re) (TLRGRADE®) (Ready-to-Use) | ALX-581-008 |
| LPS from <i>S. typhimurium</i> S-form (TLRGRADE®) (Ready-to-Use) | ALX-581-011 |
| Monophosphoryl Lipid A (Synthetic) (Ready-to-Use) | ALX-581-205 |
| Monophosphoryl Lipid A from <i>E. coli</i> , Serotype R515 (Re) (TLRGRADE®) (Ready-to-Use) | ALX-581-203 |
| Monophosphoryl Lipid A from <i>S. minnesota</i> R595 (Re) (TLRGRADE®) (Ready-to-Use) | ALX-581-202 |

FLUORESCENT DYES FOR VISUALIZING CELLULAR RESPONSES

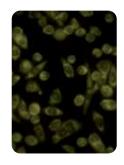
CYTO-ID® Autophagy Detection Kit 2.0 (ENZ-KIT175)

- No-transfection and quantitative approach for monitoring autophagy in live cells
- Validated under a wide range of conditions and with small molecule modulators known to influence autophagy pathways
- Stringently manufactured to control and eliminate non-specific assay artifacts



ROS-ID® Total ROS/Superoxide Detection Kit (ENZ-51010)

- Distinguishes between different reactive species such as hydrogen peroxide, peroxynitrite, and hydroxyl radicals
- High sensitivity, specificity, and accuracy for live cell studies
- Complete set of reagents, including ROS inducers and scavengers



TOOLS FOR VACCINE DEVELOPMENT

Native Bacterial and Viral Antigens

Optimized Expression Systems Deliver Superior Antigenicity

We offer a broad range of native state bacterial and viral antigens for vaccine research. Purified native protein production in the latest mammalian expression systems ensure that the resultant antigens are expressed to a high level, are in their native folded state, and possess all relevant post-translational modifications.

- High purity
- Reproducible scale-up for bulk quantities
- Lyophilized format for extended shelf-life

| BACTERIAL TOXINS AND TOXOIDS | | | | | |
|---------------------------------|-------------|--|--|--|--|
| Product Name | Product # | | | | |
| B. pertussis toxin | BML-G101 | | | | |
| B. pertussis toxin (ultra pure) | BML-G100 | | | | |
| C. difficile toxin A | BML-G140 | | | | |
| C. difficile toxin B | BML-G150 | | | | |
| C. diphtheriae toxin | BML-G135 | | | | |
| V. cholera toxin B subunit | BML-G125 | | | | |
| V. cholera toxin, (azide free) | BML-G117 | | | | |
| C. difficile toxoid A | BML-G145 | | | | |
| C. difficile toxoid B | BML-G155 | | | | |
| Tetanus toxoid | ALX-630-108 | | | | |

| RELATED PRODUCTS | | | | | | |
|---------------------------------|-------------|---------------------------------|-------------|--|--|--|
| Product Name | Product # | Product Name | Product # | | | |
| Adenovirus Type 5 hexon protein | ALX-630-120 | Dengue NS1 ST3, (recombinant) | ENZ-PRT106 | | | |
| Adenovirus type 40 antigen | ALX-630-121 | Dengue NS1 ST4, (recombinant) | ENZ-PRT107 | | | |
| CMV native antigen | ALX-630-125 | HSV-1 native antigen | ALX-630-126 | | | |
| Dengue NS1 ST1, (recombinant) | ENZ-PRT104 | HSV-2 native antigen | ALX-630-127 | | | |
| Dengue NS1 ST2, (recombinant) | ENZ-PRT105 | Yellow Fever NS1, (recombinant) | ENZ-PRT109 | | | |

Viral Agent Probes

Virus Detection Probes for use in a variety of hybridization techniques, including *in situ* hybridization (ISH) at low concentrations (50-100 ng/mL). Biotinylated probes have been shown to hybridize to homologous DNA at the same rate and to the same extent as non-biotinylated probes.

| BIO-PROBE INFECTIOUS AGENT PROBES | | | | | | |
|-----------------------------------|-----------|-------------------------------------|-----------|--|--|--|
| Product Name Product # | | Product Name | Product # | | | |
| BIO-PROBE® Adenovirus | ENZ-40834 | BIO-PROBE® Epstein-Barr Virus (EBV) | ENZ-40836 | | | |
| BIO-PROBE® BK Virus | ENZ-40848 | BIO-PROBE® Herpes Simplex | ENZ-40838 | | | |
| BIO-PROBE® Blur 8 | ENZ-40849 | BIO-PROBE® JC Virus | ENZ-40847 | | | |
| BIO-PROBE® Cytomegalovirus | ENZ-40835 | BIO-PROBE® SV 40 | ENZ-40845 | | | |

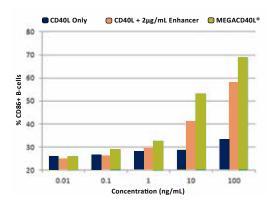
Factors for Modulating the Host Response Superior Performance from Carefully Engineered TNF Superfamily Ligands

MEGACD40L® Protein (soluble) (human), (recombinant) (ALX-522-110)

- Highly active MEGACD40L® oligomer mimics in vivo membrane-assisted CD40L aggregation and stimulation without the need for enhancers
- High purity with low endotoxin levels
- Experienced manufacturing results in consistent and dependable lot-to-lot performance
- · Optimized formulation for improved stability and enhanced activity

| RELATED PRODUCTS | | | | | |
|---|-------------|--|--|--|--|
| Product Name | Product # | | | | |
| MEGACD40L® Protein (soluble) (mouse), (recombinant) | ALX-522-120 | | | | |
| CD40L (soluble) (human), (recombinant) | ALX-522-015 | | | | |
| CD40L (soluble) (human), (recombinant) Set | ALX-850-064 | | | | |
| Fas (human):Fc (human), (recombinant) | ALX-850-016 | | | | |

Enhance Immune Activation with MEGACD40L® Protein



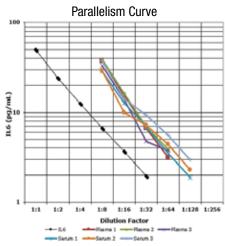
B cell lymphocyte activation by various CD40L constructs. PBMCs were treated for 48 hrs in media containing serially diluted CD40L (blue), CD40L + 2 $\mu g/mL$ Enhancer (orange), or MEGACD40L® Protein (green).

QUANTIFY BIOMARKERS OF INFLAMMATION

Detect Mediators of Acute and Chronic Inflammation

IL-6 (human), High Sensitivity ELISA Kit (ENZ-KIT178)

- Most-sensitive (0.057 pg/ mL) colorimetric kit for IL-6 quantification
- Rapid quantitative results in 3 hours
- High-throughput immunoassay measures up to 40 samples in duplicate per 96-well kit



Histamine ELISA Kit (ENZ-KIT140A)

- Highly sensitive measurement, detecting as little as 0.068 ng/mL
- Unique assay which does not require an additional acylation step
- Rapid results in 2.5 hours for up to 40 samples in duplicate

LTB₄ ELISA Kit (ADI-900-068)

- Ultrasensitive measurement of LTB4, detecting as little as 5.63 pg/ml
- Broad dynamic range suitable for a large variety of samples
- High throughput format with results in 4 hours for up to 39 samples in duplicate

| RELATED PRODUCTS | | | | | | |
|---|-------------|-------------------------|-------------|--|--|--|
| Product Name | Product # | Product Name | Product # | | | |
| Corticosterone ELISA Kit | ADI-900-097 | IFN-γ (human) ELISA Kit | ADI-900-136 | | | |
| Cortisol ELISA Kit | ADI-900-071 | IL-8 (human) ELISA Kit | ADI-900-156 | | | |
| IL-10 (human) ELISA kit | ADI-900-036 | TGF-α ELISA Kit | ENZ-KIT132 | | | |
| PGE ₂ High Sensitivity ELISA Kit | ADI-930-001 | TGF-β1 ELISA Kit | ADI-900-155 | | | |

Detect Specific Cell Populations with Antibodies for CD Markers

We offer a broad range of antibodies for use in all areas of immunology. Many are validated for use in key applications such as flow cytometry, immunohistochemistry, immunoprecipitation, and Western blotting. For more details about this program visit **www.enzolifesciences.com/antibodies**.

| CD Marker | T Cells | NK Cells | B Cells | Dendritic Cells | Granulocytes | Macrophages | Stem Cell Precursor | Endothelial Cells | Epithelial Cells |
|------------------------|---------|----------|---------|--------------------|--------------|-------------|------------------------|----------------------|---------------------|
| | | | | | | | | | |
| CD3 | • | | | | | | | | |
| CD40 | • | | • | • | • | • | • | • | • |
| CEACAM1 (CD66a) | | | | | • | | | | • |
| CEACAM6 (CD66c) | | | | | • | | | | • |
| uPAR (CD87) | • | • | | | • | • | | • | |
| CD40L (CD154) | • | | | | | | | | |
| CD163 | | | | | | • | | | |
| LAG-3 (CD223) | • | • | | | | | • | | |
| E-cadherin (CD324) | | | | | | | • | | • |
| p-glycoprotein (CD243) | | | | | | | • | | |

SCIENTISTS ENABLING HEALTHCARE™



Global Headquarters ENZO LIFE SCIENCES, INC.

81 Executive Blvd., Suite 3
Farmingdale, NY 11735
Phone: 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 Phone: +41 61 926 8989 Fax: +41 61 926 8979 info-eu@enzolifesciences.com

LOCAL EUROPEAN OFFICES

Belgium, The Netherlands & Luxembourg

Avenue Louise 65/Box 11 1050 Bruxelles Belgium Phone: +32 3 466 0420 Fax: +32 3 808 7033

Enzo Life Sciences BVBA

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 Phone: +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 Phone: +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
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
Fax: +44 1392 825910
info-uk@enzolifesciences.com

For local distributors and detailed product information visit us online: enzolifesciences.com

Simplify Your Science™ At Enzo Biochem, we focus on creating a healthier world using scientific innovation through Drug Discovery, Development, and Diagnostic products and services.

For Research Use Only. Not for use in Diagnostic Procedures. All product names, logos and brands are the property of their respective owners.



