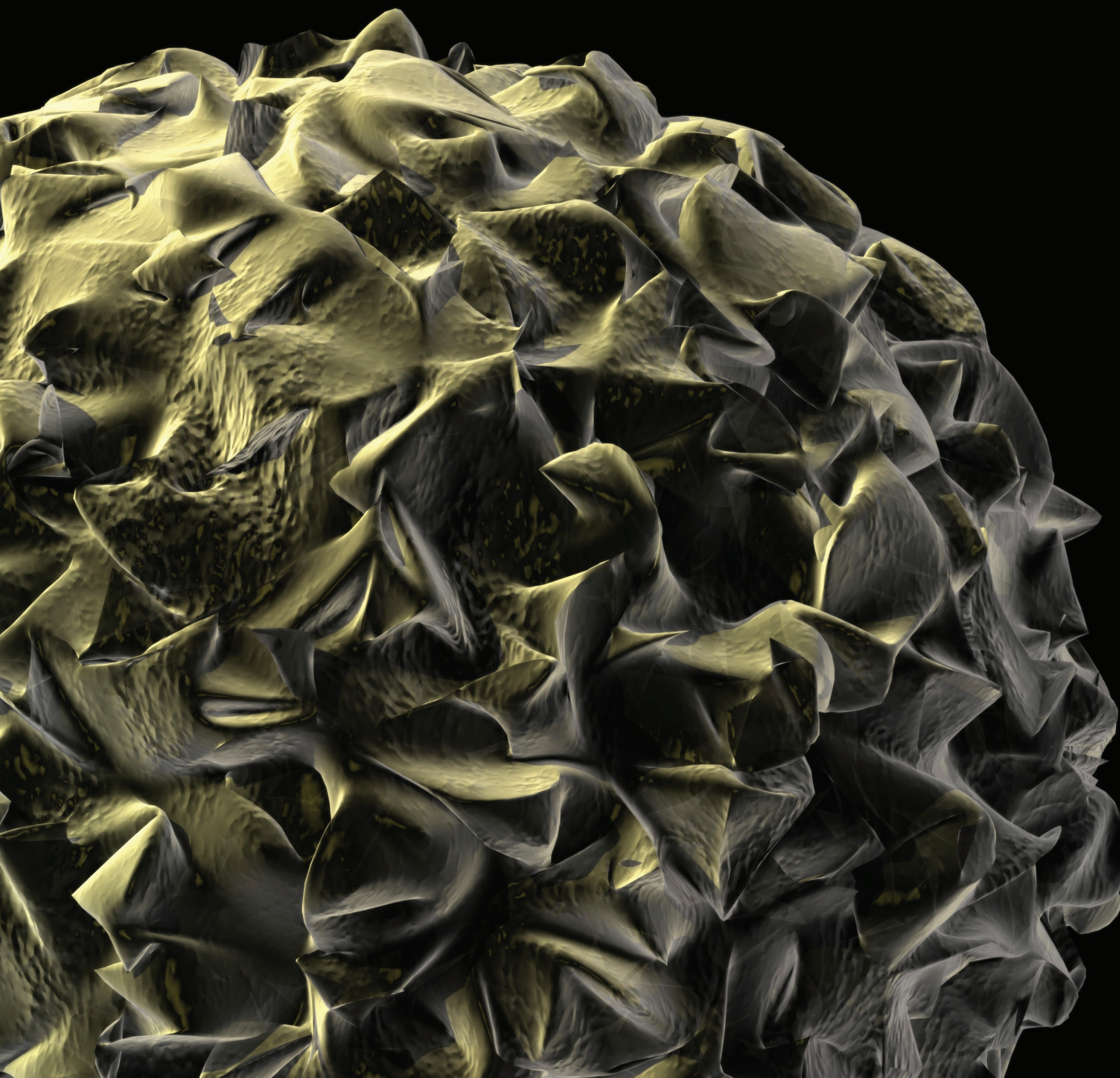




Apoptosis

Monitor Programmed Cell Death from
Membrane to Nucleus

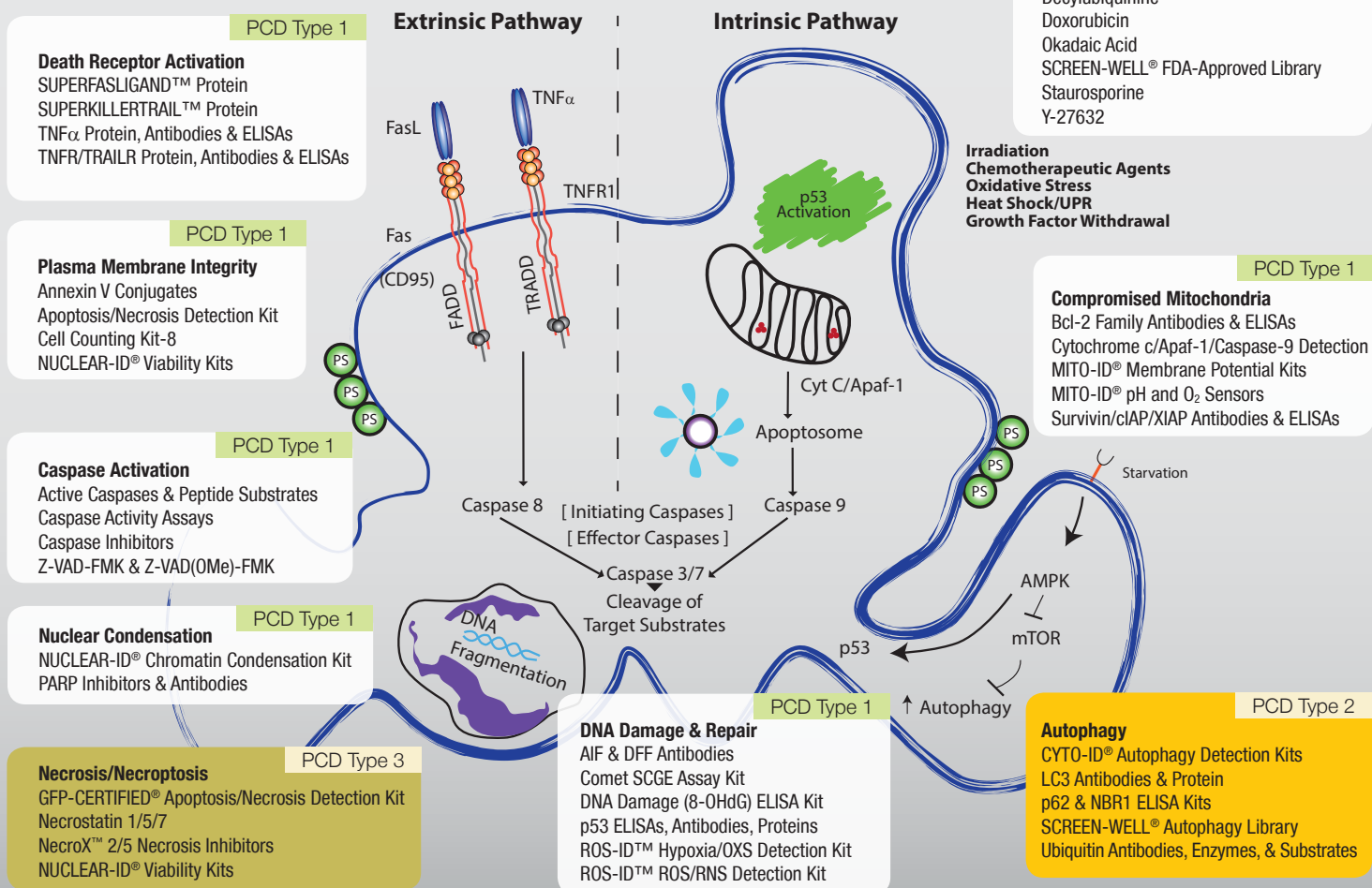


INTRINSIC. EXTRINSIC. SIMPLISTIC.

Convenient Solutions for Complete Analysis of Apoptosis and Cell Fate

Of the three major established programmed cell death (PCD) pathways, Apoptosis (PCD Type 1) is the most-well characterized, being recognized as a critical regulator of development and immunity, as well as organ and tissue homeostasis. Apoptotic cells die in a controlled fashion in response to a variety of extrinsic or intrinsic signals. Influencing cells to undergo or resist apoptosis is the aim of countless drugs for treatment of cancer, neurodegenerative, and immune-related disease. The interplay between apoptotic, autophagic, and necrotic pathways promises to yield further targets for therapeutic intervention. With over 2,000 products for the analysis of cell death, Enzo enables detection of phenotypic hallmarks of apoptotic cell death and the complete analysis of cross-talk between PCD pathways.

Apoptotic Pathways (PCD Type 1)



Live Cell Analysis

Enzo's expertise in fluorescent probe chemistry and cellular analysis combine to provide high-sensitivity fluorescent probes for characterizing cell membrane and organelle changes associated with apoptosis, including phosphatidyl serine (PS) exposure, plasma membrane integrity, mitochondrial membrane permeability, DNA damage, and chromatin condensation. Our range of unique dyes includes assays for detecting oxidative stress, ER stress, and autophagy pathways known to push cells toward an apoptotic or necrotic fate.

Immunoassays & Antibodies

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. With over 300 immunoassay and detection kits, 3,000 antibodies, and a complete set of IHC reagents, Enzo enables sensitive and specific analysis of extrinsic and intrinsic apoptosis mediators to determine the source and consequence of death pathway activation.

Modulators & Screening Assays

More than 3,000 biologically characterized small molecules, enhanced TNF ligands, and the industry's most diverse collection of SCREEN-WELL® compound libraries enables efficient induction and inhibition of apoptotic pathways. Our collection includes hundreds of caspase and protease inhibitors, FDA approved drugs, natural products, and autophagy modulators. Convenient caspase activity assays include high-purity, high-activity enzymes and cell-permeable peptide substrates and inhibitors.

DETECT HALLMARKS OF APOPTOSIS

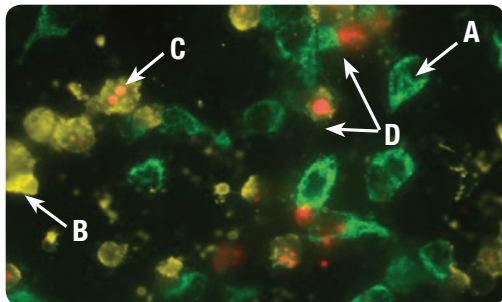
Membrane Display of Phosphatidylserine

Real-time, multiplex analysis of phosphatidylserine exposure and membrane permeability

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

- 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

Detect Cell Death in GFP-Expressing Cell Lines



GFP-CERTIFIED® Apoptosis/Necrosis Detection Kit detects four distinct cell states. Mitochondrial GFP-expressing HeLa cells were treated with 2 μ M Staurosporine for 4 hours. The Apoptosis Detection Reagent (Gold) and Necrosis Detection Reagent (Red) specifically detect cell states with clear spectral separation from mitochondria-associated GFP signal. Healthy cells (A), cells undergoing apoptosis (B), cells undergoing late-stage apoptosis (C), and necrotic cells (D).

PRODUCT NAME	PRODUCT #
Annexin V (h)(recombinant)(FITC)	ALX-209-256
Annexin V (h)(recombinant)(APC)	ALX-209-252
Annexin V (h)(recombinant)(R-PE)	ALX-209-251
Annexin V (ms)(rec.)(Alexa Fluor488)	ALX-209-257
Annexin V (h)(recombinant)(GST-tag)	ALX-201-116
Annexin V-Cy3 Apoptosis Detection Kit	ALX-850-252
Annexin V-Cy5 Apoptosis Detection Kit	ALX-850-254
Annexin V-EGFP Apoptosis Detection Kit	ALX-850-253
Annexin V-FITC Apoptosis Detection Kit	ALX-850-020
Annexin V-FITC Apoptosis Detection Kit Plus	ALX-850-255
Annexin V (h) ELISA Kit	ALX-850-049

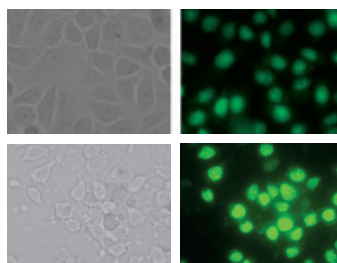
Chromatin Condensation

Sensitive, high-throughput detection of chromatin condensation

NUCLEAR-ID® Green Chromatin Condensation Kit (ENZ-51021)

- Condensation-sensitive fluorescent dye excitable with standard 488 nm laser
- Intercalating dye with superior permeability with any live cell line
- Easy no-wash, mix-and-read protocol for microplates, flow cytometry, and microscopy

Convenient alternative to Hoechst dyes that require 350nm UV laser for detection



Bright Field

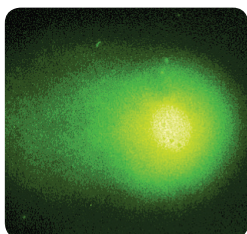
FITC Filter

Chromatin condensation as observed by fluorescence microscopy using a standard 488 nm laser. HeLa cells were treated for 4 hours with DMSO (Control) or 2 μ M Staurosporine on a slide and stained with 5 μ M NUCLEAR-ID® Green dye. The intercalating dye exhibits increased fluorescence upon chromatin condensation, a hallmark of apoptosis.

DNA Damage

Visualize DNA fragmentation in apoptotic cells

Comet SCGE Assay Kit (ADI-900-166)



- Sensitive reliable assay for detection of genotoxic agents
- Ready-to-use Comet slides allow direct application of sample without pretreatment
- Shorter assay time allows for higher throughput sample analysis

RELATED PRODUCTS

PRODUCT NAME	PRODUCT #
DNA Damage ELISA Kit	ADI-EKS-350
HSP70 High Sensitivity ELISA Kit	ADI-EKS-715
IMMUNOSET® p53/MDM2 Complex ELISA Development Set	ADI-960-070
p62 ELISA Kit	ADI-900-212
Proteasome ELISA Kit	BML-PW0575
Survivin (h) ELISA Kit	ADI-900-111

MODULATE & SCREEN CASPASE ACTIVITY

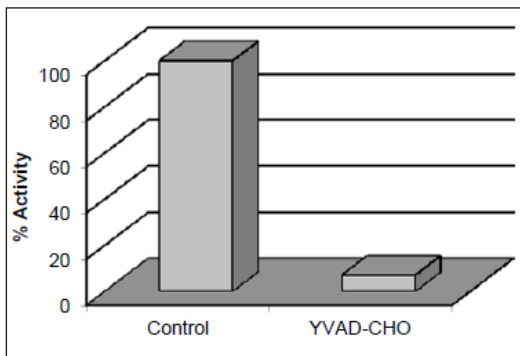
High-purity inhibitors, active enzymes, antibodies and flexible activity assays for complete analysis of caspase cascades

Caspase Activity Assays & Active Enzymes

Inhibition of Caspase-1 by Ac-YVAD-CHO

Caspase-1 Assay Kit (BML-AK701)

- Can be used to screen inhibitors of caspase-1, study enzyme regulation and kinetics, or cleave target substrates
- Flexible format allows user to choose between a colorimetric substrate or fluorogenic substrate



The enzyme was incubated with the inhibitor for 10 minutes prior to addition of substrate. 30 U/well; [YVAD-CHO]=0.1 mM; [YVAD-pNA]=200 mM; 25° C

PRODUCT NAME	PRODUCT #
Caspase Assays	
Caspase-2 Assay Kit	BML-AK702
Caspase-3 Cell Assay Kit	BML-AK703
Caspase-3 Assay Kit	BML-AK700
Caspase-8 Assay Kit	BML-AK715
Active Enzymes	
Caspase-1 (h)(recombinant)	BML-SE168
Caspase-3 (h)(recombinant)	BML-SE169
Caspase-4 (h)(recombinant)	BML-SE176
Caspase-5 (h)(recombinant)	BML-SE171
Caspase-6 (h)(recombinant)	BML-SE170
Caspase-8 (h)(recombinant)	BML-SE172

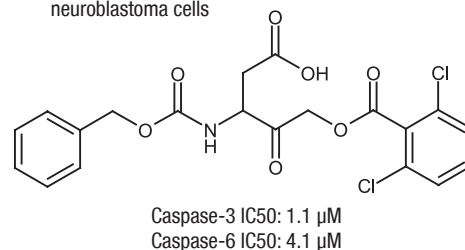
Ready-To-Use Caspase Inhibitors

High purity, peptide inhibitors for modulation of the caspase signaling cascade

PRODUCT NAME	PRODUCT #
Z-VAD-FMK [Pan-Caspase Specific Inhibitor (FMK)]	ALX-260-020
Ac-DEVD-CHO [Caspase-3/Caspase-7 Inhibitor (CHO)]	ALX-260-030
Ac-YVAD-CMK [Caspase-1 Inhibitor (CMK)]	ALX-260-028
Z-Asp-2,6-dichlorobenzoyloxymethylketone [Pan-Caspase Specific Inhibitor]	ALX-260-029
Z-DEVD-FMK (Ready-to-Use)[Caspase-3/Caspase-7 Inhibitor (FMK)]	ALX-260-141
Z-VAD-FMK (Ready-to-Use)[Pan-Caspase Specific Inhibitor (FMK)]	ALX-260-138

Z-Asp-2,6-dichlorobenzoyloxymethylketone

- Irreversible, cell permeable broad-spectrum caspase inhibitor
- Inhibits hypoxia-induced apoptosis in neuroblastoma cells



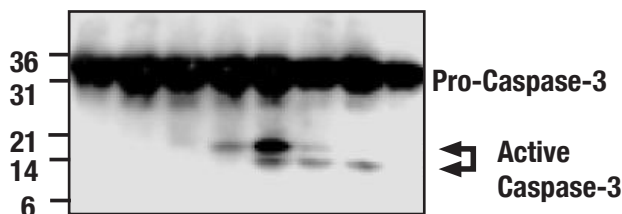
Antibody-Based Detection of Cleaved Caspases

Caspase-3, mAb (31A1067) (ALX-804-305)

- Recognizes human, mouse and rat procaspase-3 and cleaved 17 kDa catalytic subunit of caspase-3
- Validated for Western blot and immunohistochemistry applications

PRODUCT NAME	PRODUCT #
Caspase-1 pAb	ALX-210-804
Caspase-1 (ms) mAb (1H11)	ALX-804-530
Caspase-2 mAb (11B4)	ALX-804-356
Caspase-3, pAb	ADI-AAP-113
Caspase-3 mAb (31A1067)	ALX-804-305
Caspase-8 (h) mAb (12F5)	ALX-804-242
Caspase-8 mAb (5F7)	ADI-AAM-118-E

Western blot analysis for detection of Caspase-3 activation in HeLa cells



Cells were treated with 2 μM staurosporine for different time periods. Caspase-3 activation was determined by cleavage of procaspase-3, which generates 17 and 12 kDa, larger and smaller catalytic subunit, respectively.

EXTRINSIC SIGNALING THROUGH DEATH RECEPTORS

Superior performance from carefully engineered TNF Superfamily Ligands

Death Receptor Ligands

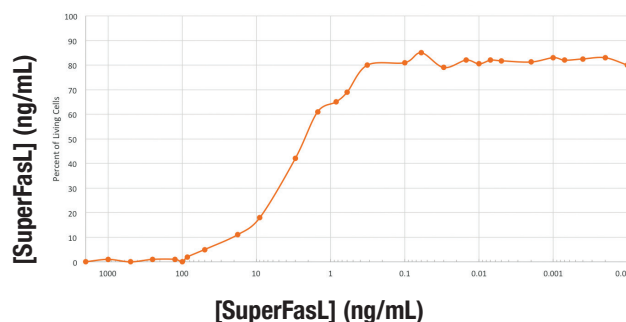
Members of the tumor necrosis factor (TNF) receptor gene superfamily bind to extrinsic ligands and transduce intracellular signals leading to the destruction of the cell. Ligands, such as FasL and TNF- α , hold potential as targets for therapeutic and diagnostic applications.

SUPERFASLIGAND™ (soluble)(human)(recombinant) (ALX-522-020)

- Features an N-Terminus linker shown to improve stability
- Enhanced immune activation compared to other recombinant ligands
- Mimics glycosylation of native human FasL

PRODUCT NAME	PRODUCT #
MEGACD40L® (soluble)(human)(recombinant)	ALX-522-110
MEGACD40L® (soluble)(mouse)(recombinant)	ALX-522-120
FasL [CD95L,CD178](soluble)(human)(recombinant)	ALX-522-001
Fc (h):TNF- α (soluble)(human)(recombinant)	ALX-522-087
KILLERTRAIL™ (soluble)(human)(recombinant)	ALX-201-073
SUPERKILLERTRAIL™ (soluble)(human)(recombinant)	ALX-201-115
SUPERKILLERTRAIL™ (soluble)(human)(recombinant)	ALX-201-130
TNF- α (human)(recombinant)(cell culture grade)	ALX-520-002
TNF- α (soluble)(human)(recombinant)	ALX-522-008
TRAIL (soluble)(human)(recombinant)(untagged)	BML-SE721
TRAIL (soluble)(ms)(recombinant)	BML-SE722

Induction of Cell Death with SUPERFASLIGAND™ Protein



Jurkat cells were incubated with the indicated concentration of SuperFasL and cell proliferation was assayed. % Viability was determined in comparison to control wells with no SuperFasL.

Oligomerization Domains in Enzo Enhanced Ligands:

Mega Ligands – use the ACRP30headless domain from human or mouse ACRP30 | **KILLER™ LIGANDS** – use a proprietary linker peptide that promotes trimerization
SUPERKILLER™ LIGANDS – use the KILLER™ linker peptide mutated to increase disulfide-mediated cross-linking

EXAMINE MITOCHONDRIAL PERTURBATIONS & INTRINSIC PATHWAY ACTIVATION

High-Throughput Assays For Mitochondrial Function

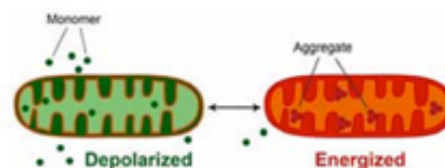
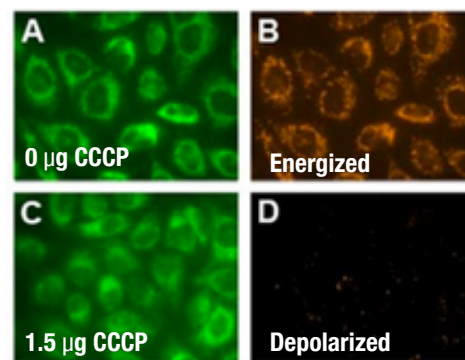
A real-time mitochondrial membrane potential assay with superior sensitivity

MITO-ID® Membrane Potential Detection Kit (ENZ-51018)

- More sensitive than JC-1
- True mix-and-read homogeneous assay for live cells
- Validated for multiple detection platforms: microscopy, flow cytometry and microplates

PRODUCT NAME	PRODUCT #
Mit-E- ψ ™ Mitochondrial Permeability Detection Kit	BML-AK116
MITO-ID® Green Detection Kit For Fluorescence Microscopy	ENZ-51022
MITO-ID® Red Detection Kit (GFP-CERTIFIED®) For Microscopy	ENZ-51007
MITO-ID® Extracellular O ₂ Sensor Kit	ENZ-51045
MITO-ID® Extracellular pH Sensor Probe	ENZ-51048

Dual-emission Dye Monitors Energetic Status



You may also be interested in:



**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

Benelux

Enzo Life Sciences BVBA
Frankrijklei 33 – Bus 31
BE-2000 Antwerpen
Belgium
Ph: +32 3 466 0420
Fax: +33 437 484 239
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: +33 437 484 239
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

Netherlands

Enzo Life Sciences BVBA
Postbus 47
NL-4940 AA Raamsdonksveer
Netherlands
Ph: +32 3 466 04 20
Fax: +33 437 484 239
info-nl@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**

Put our experience to work for you!

Our broad range of scientific expertise and industry-proven manufacturing capabilities enables us to provide a comprehensive set of apoptosis tools to save you time and money!