



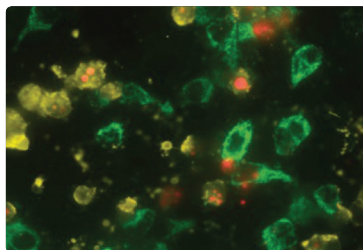
Solutions for Metal Toxicity Analysis

Innovative Metal Toxicity Assessment

In recent years, there has been increasing concern regarding metal toxicity exposure and its effect on humans. Human exposure has risen dramatically as a result of an increase of the use of harmful metals in many processes, including industrial, agricultural, domestic and technological applications. Enzo provides a comprehensive set of solutions for evaluating metal toxicity, including cell-based assays for live cell analysis, focused compound libraries for profiling organ-specific toxicity, and immunoassays for biomarker detection.

CELLESTIAL® Assays

Our CELLESTIAL® portfolio of fluorescent probes and assay kits for cellular analysis provides a complete set of tools for monitoring cell viability, proliferation, death, oxidative stress, and toxicology by flow cytometry, microscopy, and microplate platforms.



Multiplex assay that distinguishes between apoptosis and necrosis

PRODUCT NAME	PRODUCT #	PRODUCT NAME	PRODUCT #
GFP-CERTIFIED® Apoptosis/Necrosis Detection Kit	ENZ-51002	MITO-ID® Membrane Potential Cytotoxicity Kit	ENZ-51019
CYTO-ID® Autophagy Detection Kit 2.0	ENZ-KIT175	NUCLEAR-ID® Chromatin Condensation Kit	ENZ-51021
EFLUX-ID® Multidrug Resistance Kits (Gold)	ENZ-51030	PROTEOSTAT® Aggresome Detection Kit	ENZ-51035
EFLUX-ID® Multidrug Resistance Kits (Green)	ENZ-51029	ROS-ID® Total ROS Detection Kit	ENZ-51011
FLUOFORTE® Calcium Assay Kit	ENZ-51016	ROS-ID® Total ROS/Superoxide Detection Kit	ENZ-51010
LYSO-ID® Red Cytotoxicity Kit	ENZ-51015	ROS-ID® Superoxide Detection Kit	ENZ-51012

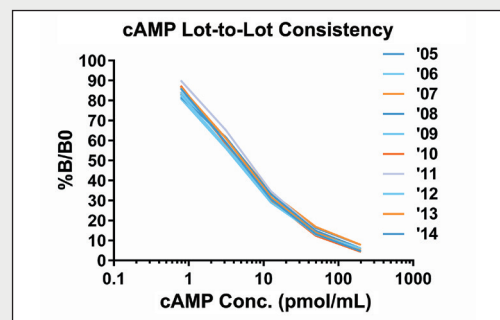
SCREEN-WELL® Compound Libraries

Enzo Life Sciences has a long and successful track record in identifying, synthesizing and commercializing known bioactives for use as research tool compounds and in assembling relevant sets of focused compounds. Our compound libraries can be used for high-throughput screening for organ-associated toxicity.

PRODUCT NAME	PRODUCT #	PRODUCT NAME	PRODUCT #
Cardiotoxicity Library	BML-2850	Myotoxicity Library	ENZ-LIB101
Hematopoietic Toxicity Library	BML-2852	Nephrotoxicity Library	ENZ-LIB100
Hepatotoxicity Library	BML-2851		

Immunoassays

We offer hundreds of ELISA kits in both immunometric and competitive assay formats. Each kit is put through rigorous testing to ensure high precision, accuracy, sensitivity, and specificity.



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

PRODUCT NAME	PRODUCT #
cAMP Complete ELISA Kit	ADI-900-163
COX-2 (human) ELISA Kit	ADI-900-094
DNA Damage ELISA Kit	ADI-EKS-350
HO-1 (human) ELISA Kit	ADI-EKS-800
HO-1 (rat) ELISA Kit	ADI-EKS-810A
KIM-1 (human) ELISA Kit	ADI-900-226
NGAL (human) ELISA Kit	BPD-KIT-036
Serotonin ELISA Kit	ADI-900-175
SMN ELISA Kit	ADI-900-209
Substance P ELISA Kit	ADI-900-018

Global Headquarters
ENZO LIFE SCIENCES, INC.
 Ph: 800.942.0430
 info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES (ELS) AG
 Ph: +41 61 926 8989
 info-eu@enzolifesciences.com

Belgium, The Netherlands & Luxembourg
& Luxembourg
 Ph: +32 3 466 0420
 info-be@enzolifesciences.com

France
 Ph: +33 472 440 655
 info-fr@enzolifesciences.com

Germany
 Ph: +49 7621 5500 526
 info-de@enzolifesciences.com

UK & Ireland
 Ph: 0845 601 1488 (UK customers)
 Ph: +44 1392 825900
 info-uk@enzolifesciences.com

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

For Research Only. Not for Use in Diagnostic Procedures.