



# Cancer Inhibitor Screening



## SCREEN-WELL® Cancer Library (ENZ-LIB102)

### Essential Standards for Cancer Inhibitor Screening

The SCREEN-WELL® Cancer Library is a collection of 275 compounds that can be used for cancer inhibitor screening and assay development. The library contains small molecules affecting mTOR, aurora kinases, tyrosine kinases, PI3K, and HDAC, as well as many structurally and mechanistically different compound classes. This library is an essential tool for cancer inhibitor screening and drug development. The compounds are dissolved in DMSO at 10 mM and aliquoted into 96-well deep-well plates at 100 µL per well.



#### PRODUCT SPECIFICATIONS

APPLICATIONS	HTS
KIT/SET CONTAINS	275 structurally and mechanistically different compounds, as well as nontoxic controls
QUANTITY	100 µL/well
CONCENTRATION	DMSO solutions (10mM)
USE/STABILITY	Stable for at least one year from the date of receipt when stored at -80° C
HANDLING	Avoid freeze/thaw cycles
SHIPPING	Shipped on dry ice
LONG TERM STORAGE	-80° C

#### RELATED PRODUCTS

APOPTOSIS/AUTOPHAGY/VIABILITY KITS	PRODUCT #
CYTO-ID® Autophagy Detection Kit 2.0	ENZ-KIT175
GFP-CERTIFIED® Apoptosis/Necrosis Detection Kit	ENZ-51002
NUCLEAR-ID® Red DNA Stain	ENZ-52406

EPIGENETIC ACTIVITY ASSAYS	PRODUCT #
CHEMILUM DE LYS® HDAC/SIRT Chemiluminescent Drug Discovery Kit	BML-AK532
FLUOR DE LYS® Green HDAC Fluorometric Activity Assay Kit	BML-AK530
FLUOR DE LYS® HDAC Fluorometric Activity Assay Kit	BML-AK500
FLUOR DE LYS® SIRT1 Fluorometric Drug Discovery Assay Kit	BML-AK555

KINASE ACTIVITY KITS	PRODUCT #
Akt Kinase Activity Kit	ADI-EKS-400A
PKA Kinase Activity Kit	ADI-EKS-390A
PKC Kinase Activity Kit	ADI-EKS-420A

SCREEN-WELL® LIBRARIES	PRODUCT #
SCREEN-WELL® FDA Approved Drug Library	BML-2843
SCREEN-WELL® Kinase Inhibitor Library	BML-2832

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  
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](http://www.enzolifesciences.com)

For Research Only. Not for Use in Diagnostic Procedures.