

# SCREEN-WELL<sup>®</sup>

## Cancer Library

### Essential standards for cancer inhibitor screening

The SCREEN-WELL<sup>®</sup> 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.

Citations: 1

[View Online »](#)

### Ordering Information

[Order Online »](#)

ENZ-LIB102-0100	100 µl/well - 1Library
-----------------	------------------------

### Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Use/Stability	Stable for at least one year from the date of receipt when stored at -80°C.
Handling	Avoid freeze/thaw cycles.
Long Term Storage	-80°C
Shipping	Dry Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Application	HTS
Concentration	DMSO solutions (10mM).
Contents	275 structurally and mechanistically different compounds, as well as nontoxic controls.
Quantity	100 µl/well

Last modified: May 29, 2024



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

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

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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