

# SCREEN-WELL®

## Kinase Inhibitor library

The SCREEN-WELL® Kinase Inhibitor Library contains 80 known kinase inhibitors of well-defined activity. Inhibitors are supplied dissolved in DMSO at 10 mM and aliquoted into deep-well plates at either 100 or 500 microliters per well. The library is an ideal tool for chemical genomics, assay development and other pharmacological applications. Includes inhibitors of these important kinases: Insulin/IGF Receptors, PI 3-Kinase, CaM Kinase II, JAK, PKA, CDK, JNK, PKC, CKI II, MAPK, RAF, EGFR, MEK, SAPK, GSK, MLCK, Src-family, IKK, PDGFR, VEGFR and many more.

Citations: 53

[View Online »](#)

### Ordering Information

[Order Online »](#)

<b>BML-2832-0100</b>	100 µl/well - 1Library
<b>BML-2832-0500</b>	500 µl/well - 1Library

### Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least one year from the date of receipt when stored at -80°C.
Long Term Storage	-80°C
Shipping	Dry Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Application	HTS
Concentration	10mM
Contents	80 compounds including inhibitors of these important kinases: BTK, CaM Kinase, CDK, CKI & II, EGFR, GSK, IKK, Insulin receptor, JAK, JNK, MAPK, MEK, MLCK, PI 3-Kinase, PDGFR, PKA, PKC, RAF, SAPK, Src-family, VEGFR, and more.
Formulation	DMSO solutions.
Quantity	100µl or 500µl per 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)