

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: 54

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-2832-0100

100 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

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

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

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

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

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