

SCREEN-WELL®

Protease Inhibitor library

The SCREEN-WELL® Protease Inhibitor Library includes 53 known protease inhibitors on a single 96-well plate. 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 includes inhibitors to a broad range of important proteases and is a convenient and cost-effective way to purchase a panel of protease inhibitors for chemical genomics, assay development, and other applications. It includes inhibitors of: Calpains, TPPII, Neutrophil Elastase, Cathepsins, DPPIV, Caspases, Granzyme B, γ -Secretase, Furin, MMPs, ACE, Kallikrein, Proteasome, Aminopeptidase B, and many more.

Citations: 7

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-2833-0100	100 μ l/well - 1Library
BML-2833-0500	500 μ l/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
-------------	-----

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)