

SCREEN-WELL[®]

Autophagy library

Structurally and mechanistically different compounds for studying the roles of pro- and anti-autophagic molecules in cells and *in vitro* applications.

The SCREEN-WELL[®] Autophagy library contains 94 compounds with defined autophagy-inducing or -inhibitory activity. Compounds are dissolved in DMSO or water at 10mM and aliquoted into deep-well plates at 100 or 500µl per well. A variety of structurally and mechanistically different compound classes are included. A useful tool for studying the roles of pro- and anti-autophagic molecules in cells as well as for use in *in vitro* applications.

Citations: 24

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-2837-0100	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.
Handling	Avoid freeze/thaw cycles.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Application	HTS
Concentration	10 mM* in DMSO***Compounds are predominantly at concentration of 10mM unless otherwise stated.**Compounds are predominantly provided in DMSO. Selected compounds are dissolved in nuclease-free water.
Contents	94 compounds with defined autophagy-inducing or -inhibitory activity. A variety of structurally and mechanistically different compound classes are included.
Quantity	100 µl/well or 500 µl/well
Technical Info / Product Notes	Application Note: <u>Cell-Based Screening of Focused Bioactive Compound Libraries: Assessing Small Molecule Modulators of the Canonical Wnt Signaling and Autophagy-Lysosome Pathways.</u>

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