SCREEN-WELL® Autophagy library

Structurally and mechanistically different compounds for studying the roles of proand 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 proand 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
BML-2837-0500	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.

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: Cell-Based Screening of Focused

Bioactive Compound Libraries: Assessing Small Molecule

Modulators of the Canonical Wnt Signaling and

Autophagy-Lysosome Pathways.

Last modified: May 29, 2024

