SCREEN-WELL[®] Wnt Pathway library

Compounds are highly relevant; many are unique

The SCREEN-WELL [®] Wnt Pathway library v 2.0 is a focused collection of 71 compounds with defined and diverse Wnt pathway activity, including activators and inhibitors of Wnts, Dishevelled, GSK-3 β and other Wnt pathway kinases, TCF/ β -catenin, DKK, LRP, Axin, Porcupine, and more. Compounds are dissolved in DMSO at 10mM or 1mM and aliquoted into deep-well plates at 100 or 500 μ l per well. A variety of structurally and mechanistically different compound classes are included. A useful tool for studying the roles of pro- and anti-Wnt pathway molecules in cells as well as for use in *in vitro* applications.

Citations: 6

View Online »

Ordering Information

Order Online »

BML-2838-0100	100 μl/well - 1Library
BML-2838-0500	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 DMSO solutions (10mM) except for Bafilomycin A which is

provided at 1mM in DMSO.

Contents 71 compounds with defined and diverse Wnt pathway

activity. A variety of structurally and mechanistically

different compound classes are included.

Quantity 100 μl/well or 500 μl/well

Last modified: May 29, 2024



eu@enzolifesciences.com