

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 μ l/well - 1Library
BML-2838-0500	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.
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



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)