

SCREEN-WELL[®] Stem Cell Library

Essential standards for stem cell regulatory and signaling pathway screening

The SCREEN-WELL[®] Stem Cell Library is a collection of 130 compounds that can be used for stem cell signaling research and drug screening. The compounds are predominantly dissolved in DMSO at 10 mM and aliquoted into 96 deep-well plates at 100 µL per well.

Ordering Information

[Order Online »](#)

ENZ-LIB103-0100	1Library
-----------------	----------

Manuals, SDS & CofA

[View Online »](#)

- Compounds can be used for diverse stem cell research to investigate cellular mechanisms that regulate self-renewal, reprogramming, and differentiation
- We provide detailed information in our comprehensive documentation for each compound



Handling & Storage

Use/Stability	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
Contents	130 structurally and mechanistically different compounds for stem cell signaling pathway research and new drug screening
Quantity	100 µl/well

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