

SCREEN-WELL®

Hepatotoxicity library

Essential standards for predictive toxicology
screening

The SCREEN-WELL® Hepatotoxicity Library is a focused collection of 238 compounds with defined and diverse hepatotoxicity, including steatosis, mitochondrial toxicity, Mallory body formation, cholestatic effects, and many more. A variety of structurally and mechanistically different compound classes are included, as well as nontoxic controls. The compounds are dissolved in DMSO at 10mM and aliquoted into deep-well plates. The library is an essential tool for predictive toxicology screening and assay development.

Citations: 6

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-2851-0100	100 µl/well - 1Library
BML-2851-0500	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).
Contents	238 compounds with defined and diverse hepatotoxicity. A variety of structurally and mechanistically different compound classes, as well as nontoxic controls, are included.
Quantity	100 µl/well or 500 µl/well

Technical Info / Product Notes

Application Note (Predictive Toxicology Application):

[Predictive High-Content/High-Throughput Assays for Hepatotoxicity Using Induced Pluripotent Stem Cell \(iPSC\)-Derived Hepatocytes.](#)
[Three Dimensional Hepatotoxicity Screening using Corning® HepatoCells, the Spheroid Microplate, and the SCREEN-WELL® Hepatotoxicity Library.](#)

Last modified: May 29, 2024

Image not found or type unknown
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