

SCREEN-WELL®

Myotoxicity library

Essential standards for predictive toxicology screening

The SCREEN-WELL® Myotoxicity Library is a focused collection of 60 compounds with defined and diverse myotoxicity, including rhabdomyolysis, cardiomyopathy, mitochondrial myopathy, vacuolar myopathy, and more. A variety of structurally and mechanistically different compound classes are included, as well as nontoxic controls. All compounds are dissolved in DMSO at 10mM unless otherwise stated in the documentation provided with the library. Selected compounds are dissolved in DI water. All compounds are provided at 100 µl volumes in a 96-well plate.

Developed in collaboration with Cytoo, a life sciences CRO focused on innovative products for the cosmetics and drug discovery markets with an emphasis in physiological cell-based assays compatible with High Content Screening, this library is an essential cost-effective tool for predictive toxicology screening and assay development.

Ordering Information

[Order Online »](#)

ENZ-LIB101-0100	100 µl/well - 1Library
-----------------	------------------------

Manuals, SDS & CofA

[View Online »](#)

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	DMSO solutions (10mM).
Contents	60 compounds with defined and diverse myotoxicity. A variety of structurally and mechanistically different compound classes, as well as nontoxic controls, are included.
Quantity	100 µl/well

Technical Info / Product Notes

Application Note:

[Higher Sensitivity of Myotoxicity Detection](#)

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)