

SCREEN- WELL® Nephrotoxicity library

Essential standards for predictive toxicology screening

The SCREEN-WELL® Nephrotoxicity Library is a focused collection of 86 compounds with defined and diverse nephrotoxicity, including analgesic nephropathy, interstitial nephritis, glomerulopathy, proximal and distal tubulopathy, acute renal failure, chronic kidney disease, and more. A variety of structurally and mechanistically different compound classes are included, as well as nontoxic controls. Most of the 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. This library is an essential cost-effective tool for predictive toxicology screening and assay development.

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

ENZ-LIB100-0100	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	Solvent and concentration are specified in the documentation provided with the library.
Contents	86 compounds with defined and diverse nephrotoxicity. A variety of structurally and mechanistically different compound classes, as well as nontoxic controls, are included.
Quantity	100 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)