# SCREENWELL® 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, glomerularopathy, 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 µI 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



info-

eu@enzolifesciences.com