

# SCREEN-WELL®

## Hematopoietic Toxicity library

### Essential standards for predictive toxicology screening

The SCREEN-WELL® Hematopoietic Toxicity Library is a focused collection of 115 compounds with defined and diverse hematopoietic toxicity, including myelosuppression, neutropenia, leukopenia, anemia, 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.

### Ordering Information

[Order Online »](#)

BML-2852-0100	100 µl/well - 1Library
---------------	------------------------

### Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

<b>Use/Stability</b>	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.
<b>Handling</b>	Avoid freeze/thaw cycles.
<b>Long Term Storage</b>	-80°C
<b>Shipping</b>	Dry Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

<b>Application</b>	HTS
<b>Concentration</b>	DMSO solutions (10mM).
<b>Contents</b>	115 compounds with defined and diverse hematopoietic toxicity. A variety of structurally and mechanistically different compound classes, as well as nontoxic controls, are included.
<b>Quantity</b>	100 µl/well or 500 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)