



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 |
| BML-2852-0500 | 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 | 115 compounds with defined and diverse hematopoietic toxicity. A variety of structurally and mechanistically different compound classes, as well as nontoxic controls, are included. |
| Quantity | 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)