

SCREEN-WELL®

REDOX library

The SCREEN-WELL® REDOX library contains 83 compounds with defined prooxidant or antioxidant activity. Compounds are dissolved in DMSO at 10mM and aliquoted into deep-well plates at 100 or 500µl per well. A variety of structurally and mechanistically different compound classes are included such as hydroperoxides, polyphenolics, metal chelators, thiols, thiol traps, radical scavengers, lazarooids and glutathione modulators as well as small molecule enzyme mimetics for SOD and glutathione peroxidase. The library is a useful tool for studying the roles of pro- and antioxidant molecules in cells as well as for use in *in vitro* applications.

Citations: 17

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-2835-0100	100 µl/well - 1Library
BML-2835-0500	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.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Application	HTS
-------------	-----

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)