

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, lazaroids 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

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

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com