

SCREEN-WELL®

Nuclear Receptor ligand library

The SCREEN-WELL® Nuclear Receptor Ligand Library contains 74 compounds with defined, putative and potential activity at nuclear receptors. Receptor agonists and antagonists are included. The library consists of a chemically diverse group of compounds which are supplied dissolved in DMSO at 10 mM (1mM for 5 compounds) and aliquoted into deep-well, 96-well plates at 500 µl or 100 µl per well. The library is an ideal tool for chemical genomics, receptor de-orphaning and other routine pharmacological applications. The library includes ligands for the following classes of nuclear receptors: AHR, CAR, ER, FXR, LXR, PPAR, PXR, RAR, RXR, VDR, and other steroid receptors.

Citations: 16

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-2802-0100	100 µl/well - 1Library
BML-2802-0500	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

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