

SCREEN-WELL®

Endocannabinoid library

The SCREEN-WELL® Endocannabinoid Library contains 58 compounds with defined, putative, potential or speculative activity at cannabinoid (CB) and vanilloid (VR) receptors. The library is an array of 10 different fatty acids and 6 different polar head groups. Ligands are supplied dissolved in biocompatible solvents at 1 mM and aliquoted into deep-well plates at 100 µl per well. The library is a rich source of ligands for receptor de-orphaning. Includes the following classes of ligands: Amides, Ethanolamides, Lip amino acids, Acyl-GABAs, Acyl-dopamines.

Citations: 4

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-2801-0500	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