

SCREEN-WELL®

Bioactive lipid library

The SCREEN-WELL® Bioactive Lipid Library contains 190 bioactive lipids in a 96 well format. Ideal for screening or identifying recombinant orphan G protein-coupled receptors, target validation, secondary screening, validating new assays and for routine pharmacological applications.

Citations: 20

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-2800-0100	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
Concentration	DMSO solutions (1.0 or 0.1mM, depending on compound).
Contents	190 compounds including: Agonists & Antagonists, Endocannabinoids, Farnesyl/geranylgeranyl derivatives, HETEs deHETEs, hepoxilins, Polyunsaturated fatty acids, Leukotrienes, lipoxins, LPA & phosphatidic acids, Octadecanoids, PAFs, Prostaglandins & thromboxanes, Retinoids, vitamin D metabolites, Sphingolipids.
Quantity	100 or 500µl per well.

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