

# 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-2800-0500	500 µl/well - 1Library
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](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)