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, Lipoamino acids, Acyl-GABAs, Acyl-dopamines.

Citations: 4

View Online »

Ordering Information

Order Online »

BML-2801-0500	500 μl/well - 1Library
BML-2801-0100	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



eu@enzolifesciences.com