

SCREEN-WELL® Fatty Acid library

The SCREEN-WELL® Fatty Acid Library contains 64 fatty acids of varying saturation and chain length. Many free fatty acids and their metabolites serve as precursors to important signaling systems (e.g. prostaglandins), act directly as receptor ligands or enzyme inhibitors, and have pro- or anti-atherosclerotic and anti-hypertensive properties through mechanisms that are still being characterized. Therefore, a diverse set of fatty acids can be a rich source of compounds for receptor de-orphaning, inhibitor screening, and high content screening. Fatty acids are supplied dissolved in DMSO at 10 mM and aliquoted into deep-well plates at either 100 or 500 µl per well. This set contains both fatty acids with defined activities and fatty acids with putative biological activities. Library includes: Unsaturated & saturated fatty acids, Arachidonic acid, C10-24 chain length Linolenic acid, E and Z isomers, γ-Linolenic acid and many more.

Citations: 11

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-2803-0100

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	10mM in DMSO

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