

Screen-Well™ Fatty Acid Library

CATALOG NO.: 2803

Version: 1.0

Lot #: SAMPLE

DESCRIPTION: The Screen-Well™ Fatty Acid Library contains 68 fatty acids of varying saturation and chain length. Many 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 or anti-hypertensive properties through mechanisms that are still being characterized. This library contains both fatty acids with defined activities and fatty acids with putative biological activities. Useful for receptor de-orphaning, inhibitor screening and high content screening.

CONTENTS: 68 compounds including

- Unsaturated and saturated fatty acids
- C10-24 chain length
- E and Z isomers
- Arachidonic acid
- Linolenic Acid
- γ-Linolenic acid
- ...and more

The complete list of compounds is available within the documentation included with the library or by request.

FORMAT: 10 mM DMSO solutions. 1 x 96-well plate. 100 or 500µl per well.

SUPPORTING INFORMATION: Documentation containing a complete list of compounds and supporting information is supplied with the library or by request at compoundlibraries@enzolifesciences.com. Documentation includes physical information, plate positions, and a structure database (SD) file for integration into your compound database.

STORAGE: -80°C. Thaw the plates in a 37°C bath.

FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.