N-Arachidonoyl-L-phenylalanine

Putative endogenous member of a growing family of lipo-amino acids (N-acyl amino acids). Product is subject to oxidation.

Citations: 1

View Online »

Ordering Information

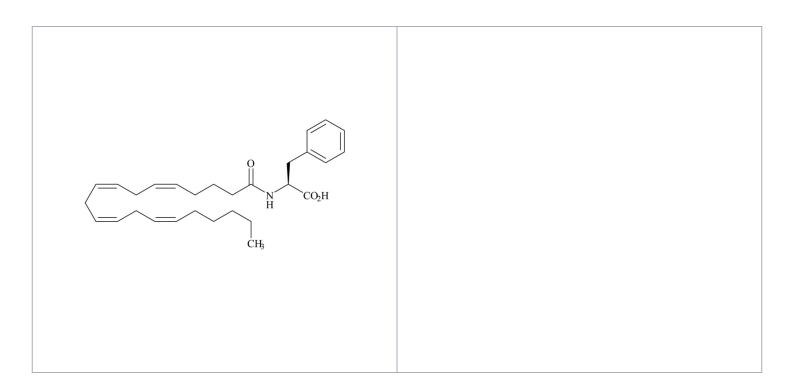
Order Online »

BML-FA040-0025

25mg

Manuals, SDS & CofA

View Online »



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended. Stable for at least

1 year after receipt when stored, as supplied, at -80°C and under an inert atmosphere.

Store solutions at -80°C for up to one month.

Handling Keep under inert gas.

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Appearance Pale yellow viscous oil.

Formula $C_{29}H_{41}NO_3$

MW 451.6

Purity ≥98% (TLC)

Soluble in DMSO (>25mg/ml) or 100% ethanol

(>25mg/ml).

Last modified: May 29, 2024



eu@enzolifesciences.com