

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 »](#)

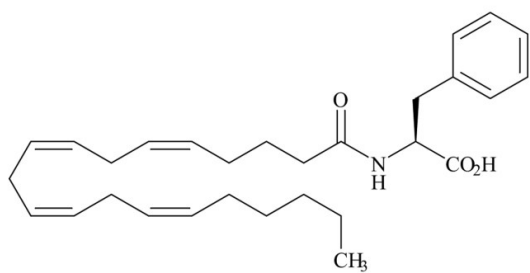
Ordering Information

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

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)