

Linolenic acid (18:3, n-3)

PPAR ligand

Useful for inducing adipocyte differentiation and/or activating PPARs in cultured cells, e.g. 200 and 250µM linolenic acid for, respectively, stimulation of FA transporter transcription in FAAR (NUC I) transfected 3T3-C2 cells and differentiation of Ob1771 preadipocytes. The product is not sterile.

Citations: 1

[View Online »](#)

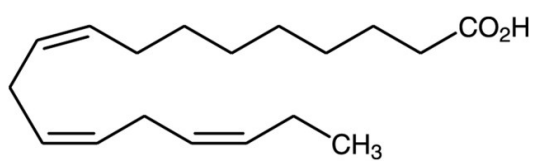
Ordering Information

[Order Online »](#)

BML-FA005-1000	1g
----------------	----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from air.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Octadeca-9Z,12Z,15Z-trienoic acid
Appearance	Near-colorless oil.
CAS	463-40-1
Couple Target	PPAR
Couple Type	Ligand
Formula	$C_{18}H_{30}O_2$
MW	278.4
Purity	≥99% (GC)
Solubility	Soluble in ethanol, hexane and methylene chloride.

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-be@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-fr@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)