

Eicosapentaenoyl ethanolamide

Has been shown to inhibit dietary-restriction-induced lifespan extension in wild-type and TOR pathway mutant nematodes. This data indicates that reduced N-acylethanolamide signaling mediates some of the effect of dietary-restriction-induced lifespan extension and that EPEA, specifically, acts as a metabolic signal that couples nutrient availability with growth and lifespan. EPEA has also been shown to have anti-proliferative effects and anti-inflammatory properties. Compound is subject to oxidation. Protect from air.

Citations: 3

[View Online »](#)

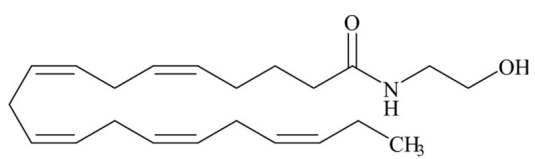
Ordering Information

[Order Online »](#)

BML-FA047-0025	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 one year after receipt when stored at -20°C. Solutions can be stored at -20°C for up to 3 months.
Handling	Protect from air.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	EPEA
Appearance	Pale amber oil.
CAS	109001-03-8
Formula	$C_{22}H_{35}NO_2$
MW	345.5
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (20mg/ml) or in 100% ethanol (25mg/ml).
Source	Synthetic.



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)