

# 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 »](#)

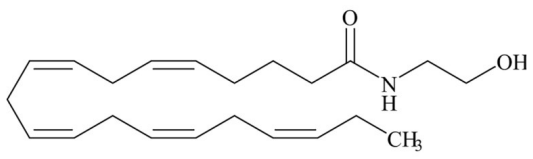
## Ordering Information

[Order Online »](#)

BML-FA047-0025	25mg
----------------	------

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

<b>Use/Stability</b>	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.
<b>Handling</b>	Protect from air.
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

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

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)