

Eicosapentaenoic acid

Inhibits PGE₂ formation

Beneficial effect in cardiovascular diseases and inflammation. Inhibits PGE₂ formation from arachidonic acid in 3T3 fibroblasts.

Citations: 4

[View Online »](#)

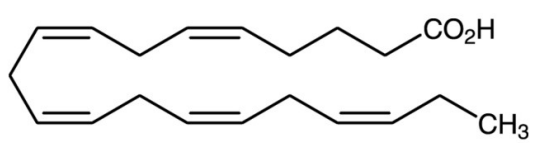
Ordering Information

[Order Online »](#)

BML-FA001-0100	100mg
BML-FA001-1000	1g

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 -20°C. After opening, store under an inert atmosphere (e.g. N2 or argon) at -80°C. Stock solutions are stable for up to 2 months at -80°C under an inert atmosphere.
Handling	Protect from air.
Long Term Storage	-20°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Eicosa-5Z,8Z,11Z,14Z,17Z-pentaenoic acid (20:5, n-3), Icosapent
Appearance	Colorless oil.
CAS	10417-94-4
Couple Target	Prostaglandin E2
Couple Type	Inhibitor
Formula	$C_{20}H_{30}O_2$
MW	302.5
Purity	≥98% (TLC)
Solubility	Soluble in 100% ethanol.

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-
fr@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)