

Arachidonic acid

(20:4, n-6)

Eicosanoid precursor and second messenger

Arachidonic acid is an essential fatty acid and a precursor to a large family of eicosanoids. Arachidonic acid acts as a second messenger independent of metabolism and a retrograde messenger in long-term potentiation in the nervous system. Arachidonic acid activates several protein kinase C (PKC) isotypes and inhibits ras-GAP.

Citations: 23

[View Online »](#)

Ordering Information

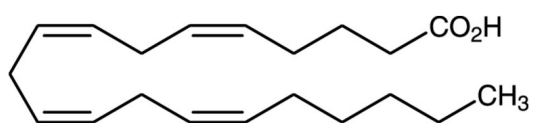
[Order Online »](#)

BML-FA003-0100	100mg
BML-FA003-1000	1g

Manuals, SDS & CofA

[View Online »](#)

- Essential precursor for leukotrienes, prostaglandins, and thromboxanes
- Used in cancer, cardiovascular, cell signaling, immunology, and metabolism studies amongst others



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for 12 months after receipt when stored at -20°C. Keep aqueous solutions on ice and use within 12 hours.
Handling	Protect from light and oxygen.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	5Z,8Z,11Z,14Z-Eicosatetraenoic acid
Appearance	Colorless oil.
CAS	506-32-1
Couple Target	PKC, Ras-GAP
Couple Type	Activator, Inhibitor
Formula	$C_{20}H_{32}O_2$
MI	14: 765
MW	304.5
Purity	≥99% (GLC, TLC)
RTECS	CE6675000
Solubility	Soluble in DMF (100mg/ml), DMSO (100mg/ml), or 100% ethanol (50mg/ml). Sparingly soluble in basic aqueous buffers (1.7mg/ml). To prevent oxidation, the solvent should be purged with an inert gas.



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)

Last modified: May 29, 2024