

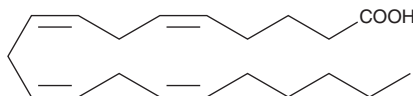
Product Information



Arachidonic Acid Lipid Maps MS Standard

Catalog No. 10007268

CAS Registry No.: 506-32-1
Formal Name: 5Z,8Z,11Z,14Z-Eicosatetraenoic acid
MF: C₂₀H₃₂O₂
FW: 304.5
Purity: ≥98%
Stability: ≥1 year at -20°C
Supplied as: A solution in ethanol



Laboratory Procedures

For long term storage, we suggest that arachidonic acid be stored as supplied at -20°C. It will be stable for at least one year.

Arachidonic acid is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as DMSO or dimethyl formamide can be used. To prevent oxidation of arachidonic acid, the solvent should be purged with an inert gas. The solubility of arachidonic acid in these solvents is at least 100 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. If an organic solvent-free aqueous solution of arachidonic acid is needed, it can be prepared by evaporating the ethanol and directly dissolving the neat oil in basic buffers. Arachidonic acid is sparingly soluble in neutral buffers. Store aqueous solutions of arachidonic acid on ice and use within 12 hours of preparation. Although the aqueous solutions of arachidonic acid may be stable for more than 12 hours, we strongly recommend using a fresh preparation each day.

Arachidonic acid is an essential fatty acid and a precursor for all prostaglandins, thromboxanes, and leukotrienes. Virtually all cellular arachidonic acid is esterified in membrane phospholipids where its presence is tightly regulated through multiple interconnected pathways.¹ Free arachidonic acid is a transient, critical substrate for the biosynthesis of eicosanoid second messengers. Receptor-stimulated release, metabolism, and re-uptake of free arachidonate are all important aspects of cell signaling and inflammation.²

References

1. Nixon, A.B., Greene, D.G., and Wykle, R.L. Comparison of acceptor and donor substrates in the CoA-independent transacylase reaction in human neutrophils. *Biochim. Biophys. Acta* **1300**, 187-196 (1996).
2. Burgoyne, R.D. and Morgan, A. The control of free arachidonic acid levels. *Trends Biochem. Sci.* **15**, 365-366 (1990).

Related Products

Arachidonic Acid - Cat. No. 90010 • Arachidonic Acid methyl ester - Cat. No. 90014 • Arachidonic Acid-d8 - Cat. No. 390010 • *trans*-4,5-epoxy-2(E)-Decenal - Cat. No. 10004257 • DDA - Cat. No. 10005432 • Methyl Pentacosanoate - Cat. No. 10006451 • Dihomo- γ -Linolenic Acid methyl ester - Cat. No. 10006580 • Arachidonic Acid (sodium salt) - Cat. No. 10006607 • Arachidonic Acid Quant-PAK - Cat. No. 10006835 • Arachidonic Acid-biotinamide - Cat. No. 10007466 • 20-carboxy Arachidonic Acid - Cat. No. 10007912 • Arachidonic Acid ethyl ester - Cat. No. 10008200 • Arachidonic Acid Leelamide - Cat. No. 10008617 • 15-Lipoxygenase-2 (human recombinant) - Cat. No. 10011263 • 15-Lipoxygenase-2 Western Ready Control - Cat. No. 10011500

Cayman Chemical

Mailing address
1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone
(800) 364-9897
(734) 971-3335

Fax
(734) 971-3640

E-Mail
custserv@caymanchem.com

Web
www.caymanchem.com

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY; NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent under separate cover to the MSDS supervisor at your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery**.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees. Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy located on our website and in our catalog**.

Copyright Cayman Chemical Company, 07/06/2009