

Product Information

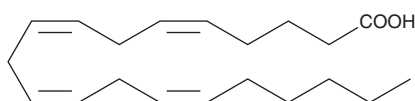


Arachidonic Acid Quant-PAK

Catalog No. 10006835

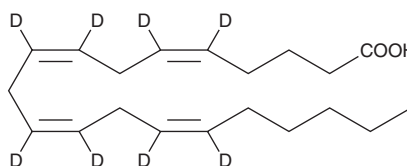
Arachidonic Acid

CAS Registry No.: 506-32-1
Formal Name: 5Z,8Z,11Z,14Z-eicosatetraenoic acid
MF: C₂₀H₃₂O₂
FW: 304.5
Purity: ≥99%
Stability: ≥1 year at -20°C
Supplied as: A solution in ethanol



Arachidonic Acid-d₈

CAS Registry No.: 69254-37-1
Formal Name: 5Z,8Z,11Z,14Z-eicosatetraenoic-5,6,8,9,11,12,14,15-d₈ acid
MF: C₂₀H₂₄D₈O₂
FW: 312.5
Chemical Purity: ≥98%
Deuterium Incorporation: ≤1% d₀
Stability: ≥1 year at -20°C
Supplied as: A solution in methyl acetate



This arachidonic acid Quant-PAK contains 50 µg of arachidonic acid-d₈ and 2-4 mg of arachidonic acid (please see the vial for exact amount and concentration). For long term storage, we suggest that arachidonic acid and arachidonic acid-d₈ be stored as supplied at -20°C. They should 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 and dimethyl formamide purged with an inert gas can be used. The solubility of arachidonic acid in these solvents is approximately 100 mg/ml.

Arachidonic acid-d₈ is supplied as a solution in methyl acetate. To change the solvent, simply evaporate the methyl acetate under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide (DMF) purged with an inert gas can be used. Arachidonic acid-d₈ is miscible in ethanol and has a solubility of approximately 100 mg/ml in DMSO and DMF.

Arachidonic acid-d₈ contains eight deuterium atoms at the 5, 6, 8, 9, 11, 12, 14, and 15 positions. It is intended for use as an internal standard for the quantification of arachidonic acid-d₈ by GC- or LC-mass spectrometry. The accuracy of the sample weight in the arachidonic acid-d₈ vial is between 5% over and 2% under the weight indicated on the vial. For better precision we have provided a precisely weighed unlabeled arachidonic acid, with the precise weight (2-4 mg) indicated on the vial. Using this vial the deuterated standard can be quantified by constructing a standard curve of peak intensity ratios (deuterated *versus* unlabeled).

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 signalling and inflammation.²

References

1. Nixon, A.B., Greene, D.G., and Wykle, R.L. *Biochim. Biophys. Acta* **1300**, 187-196 (1996).
2. Burgoyne, R.D. and Morgan, A. *Trends Biochem. Sci.* **15**, 365-366 (1990).

Related Products

Arachidonic Acid - Cat. No. 90010 • Arachidonic Acid-d₈ - Cat. No. 390010 • Docosahexaenoic Acid Quant-PAK - Cat. No. 10006829 • ω-3 Arachidonic Acid Quant-PAK - Cat. No. 10006833 • Linoleic Acid Quant-PAK - Cat. No. 10006834

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 *via* email to 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, 02/11/2010

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