

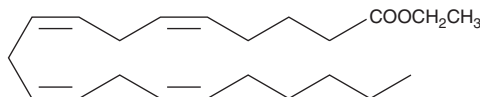
PRODUCT INFORMATION



Arachidonic Acid ethyl ester

Item No. 10008200

CAS Registry No.: 1808-26-0
Formal Name: 5Z,8Z,11Z,14Z-eicosatetraenoic acid, ethyl ester
Synonym: Ethyl Arachidonate
MF: $C_{22}H_{36}O_2$
FW: 332.5
Purity: $\geq 98\%$
Supplied as: A solution in ethanol
Storage: $-20^{\circ}C$
Stability: ≥ 1 year



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Arachidonic acid ethyl ester 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 ethyl ester in these solvents is approximately 100 mg/ml.

Description

Arachidonic acid is the keystone essential fatty acid at the origin of the arachidonic acid cascade. It is converted by cyclooxygenase, lipoxygenase, and epoxygenase enzymes into more than one hundred fifty different potent primary autacoid metabolites in species ranging from fungi to plants to mammals. Arachidonic acid is stored in tissue phospholipids in esterified form, where it comprises a small but critically controlled percentage of the polyunsaturated fatty acid pool.¹ Arachidonic acid ethyl ester is a more lipophilic form of arachidonic acid that can be incorporated into dietary regimens or fed to cultured cells as a source of exogenous arachidonate. It is one of the fatty acid ethyl esters that increase cytosolic Ca^{2+} concentration leading to pancreatic acinar cell injury due to excessive consumption of ethanol.² Whereas arachidonic acid inhibits dopamine uptake, the ethyl esterified version does not retain this property.³

References

1. Holman, R.T. Control of polyunsaturated acids in tissue lipids. *J. Am. Coll. Nutr.* **5(2)**, 183-211 (1986).
2. Chen, N., Appell, M., Berfield, J.L., *et al.* Inhibition by arachidonic acid and other fatty acids of dopamine uptake at the human dopamine transporter. *Eur. J. Pharmacol.* **478(2-3)**, 89-95 (2003).
3. Criddle, D.N., Raraty, M.G.T., Neoptolemos, J.P., *et al.* Ethanol toxicity in pancreatic acinar cells: Mediation by nonoxidative fatty acid metabolites. *Proc. Natl. Acad. Sci. USA* **101(29)**, 10738-10743 (2004).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/15/2021

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM