

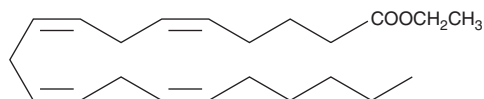
PRODUCT INFORMATION



Arachidonic Acid ethyl ester

Item No. 10008200

CAS Registry No.: 1808-26-0
Formal Name: 5Z,8Z,11Z,14Z-eicosatetraenoic acid, ethyl ester
Synonyms: C20:4 (cis-5,8,11,14) ethyl ester, Ethyl Arachidonate, Ethyl (cis-5,8,11,14)-eicosatetraenoate, SFE 22:4
MF: C₂₂H₃₆O₂
FW: 332.5
Purity: ≥98%
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥2 years



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 ethyl ester is an esterified form of arachidonic acid (Item Nos. 90010 | 90010.1 | 10006607). It is commonly used as a source of exogenous arachidonic acid *in vivo*.¹ Arachidonic acid ethyl ester has been found in the meconium of newborns suspected of fetal alcohol exposure.²

References

1. Naoe, S., Tsugawa, H., Takahashi, M., *et al.* Characterization of lipid profiles after dietary intake of polyunsaturated fatty acids using integrated untargeted and targeted lipidomics. *Metabolites* **9**(10), 241 (2019).
2. Moore, C., Jones, J., Lewis, D., *et al.* Prevalence of fatty acid ethyl esters in meconium specimens. *Clin. Chem.* **49**(1), 133-136 (2003).

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, 03/27/2024

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