

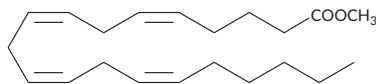
# PRODUCT INFORMATION



## Arachidonic Acid methyl ester

Item No. 90014

**CAS Registry No.:** 2566-89-4  
**Formal Name:** 5Z,8Z,11Z,14Z-eicosatetraenoic acid, methyl ester  
**Synonyms:** C20:4 (cis-5,8,11,14) methyl ester, Methyl Arachidonate, Methyl (cis-5,8,11,14)-eicosatetraenoate, SFE 21:4  
**MF:** C<sub>21</sub>H<sub>34</sub>O<sub>2</sub>  
**FW:** 318.5  
**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 methyl ester is supplied as a solution in ethanol. To change the solvent, simply evaporate the arachidonic acid methyl ester under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide purged with an inert gas can be used. Arachidonic acid methyl ester is miscible in these solvents.

### Description

Arachidonic acid methyl ester is an esterified form of arachidonic acid (Item Nos. 90010 | 90010.1 | 10006607). It is commonly used as a reference standard for the quantification of arachidonic acid in biological samples and as a source of exogenous arachidonic acid in cells and *in vivo*.<sup>1-4</sup>

### References

1. Leichsenring, M., Sütterlin, N., Less, S., *et al.* Polyunsaturated fatty acids in erythrocyte and plasma lipids of children with severe protein-energy malnutrition. *Acta Paediatr.* **84(5)**, 516-520 (1995).
2. Burdge, G.C. and Postle, A.D. Phospholipid molecular species composition of developing fetal guinea pig brain. *Lipids* **30(8)**, 719-724 (1995).
3. Martins, A.P., Yokoya, N.S., and Colepicolo, P. Comparison of extraction and transesterification methods on the determination of the fatty acid contents of three Brazilian seaweed species. *Braz. J. Pharmacol.* **22(4)**, 854-860 (2012).
4. Laborit, H. and Valette, N. The action of arachidonic acid on experimental hypertension in the rat. *Chem. Biol. Interact.* **10(4)**, 239-246 (1975).

#### 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, 02/14/2024

#### 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