

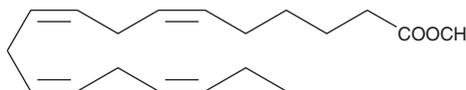
PRODUCT INFORMATION



Stearidonic Acid methyl ester

Item No. 10005000

CAS Registry No.: 73097-00-4
Formal Name: 6Z,9Z,12Z,15Z-octadecatetraenoic acid, methyl ester
Synonyms: Morocitic Acid methyl ester, SFE 19:4
MF: C₁₉H₃₀O₂
FW: 290.4
Purity: ≥98%
Supplied as: A solution in methanol
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

Stearidonic acid methyl ester is supplied as a solution in methanol. To change the solvent, simply evaporate the methanol 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. The solubility of stearidonic acid methyl ester in these solvents is approximately 50 mg/ml.

Stearidonic acid methyl ester is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the methanolic solution of stearidonic acid methyl ester should be diluted with the aqueous buffer of choice. The solubility of stearidonic acid methyl ester in PBS (pH 7.2) is approximately 0.15 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Stearidonic acid (18:4 ω -3) is a rare polyunsaturated fatty acid of the ω -3 series. In normal humans, it represents less than 0.25% of serum phospholipid fatty acids.¹ It is present in certain natural oils such as echium and black currant seed, and to the extent that these oils are incorporated into nutraceuticals, stearidonic acid can become a significant polyunsaturate in the human diet. Stearidonic acid methyl ester is an ester version of the free acid that is less water soluble, but more amenable for the formulation of stearidonate-containing diets and dietary supplements.

References

1. Holman, R.T. Control of polyunsaturated acids in tissue lipids. *J. Am. Coll. Cardiol.* **5**, 183-211 (1986).

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, 03/13/2024

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