

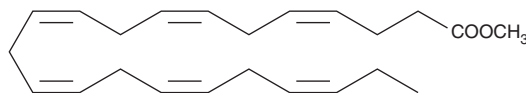
PRODUCT INFORMATION



Docosahexaenoic Acid methyl ester

Item No. 10006865

CAS Registry No.: 2566-90-7
Formal Name: 4Z,7Z,10Z,13Z,16Z,19Z-docosahexaenoic acid, methyl ester
Synonyms: Cervonic Acid methyl ester, C22:6 (cis-4,7,10,13,16,19) methyl ester, DHA methyl ester, all-cis-4,7,10,13,16,19-DHA methyl ester, Methyl all-cis-4,7,10,13,16,19-Docosahexaenoate
MF: C₂₃H₃₄O₂
FW: 342.5
Purity: ≥98%
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥1 year



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

Laboratory Procedures

Docosahexaenoic acid methyl ester (DHA methyl 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 DHA methyl ester in these solvents is approximately 100 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. If an organic solvent-free solution of DHA methyl ester is needed, it can be prepared by evaporating the ethanol and directly dissolving the neat oil in aqueous buffers. The solubility of DHA 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

Docosahexaenoic acid (DHA) methyl ester is an esterified form of DHA (Item No. 90310). It has been used as a reference standard in the quantification of fatty acids in microalgal and fish oils.¹

Reference

1. Armenta, R.E., Scott, S.D., Burja, A.M., *et al.* Optimization of fatty acid determination in selected fish and microalgal oils. *Chromatographia* **70**(3), 629–636 (2009).

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, 07/08/2020

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