

PRODUCT INFORMATION



Docosapentaenoic Acid methyl ester

Item No. 9001870

CAS Registry No.: 108698-02-8
Formal Name: (7Z,10Z,13Z,16Z,19Z)-7,10,13,16,19-docosapentaenoic acid, methyl ester

Synonyms: C22:5 (cis-7,10,13,16,19) methyl ester, all-cis-7,10,13,16,19-DPA methyl ester, all-Z-7,10,13,16,19-DPA methyl ester, DPA methyl ester, Methyl all-cis-7,10,13,16,19-Docosapentaenoate, n-3 DPA methyl ester, SFE 23:5

MF: C₂₃H₃₆O₂

FW: 344.5

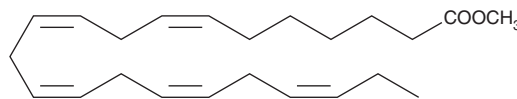
Purity: ≥98%

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

Docosapentaenoic acid 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, dimethyl formamide (DMF), and 0.1 M Na₂CO₃ purged with an inert gas can be used. The solubility of docosapentaenoic acid methyl ester in DMSO and DMF is approximately 100 mg/ml and approximately 1 mg/ml in 0.1 M Na₂CO₃.

Description

Docosapentaenoic acid methyl ester is an esterified form of docosapentaenoic acid (Item Nos. 90165 | 21907). 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, 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