

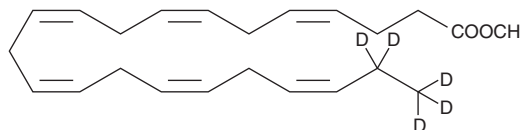
PRODUCT INFORMATION



Docosahexaenoic Acid-d₅ methyl ester

Item No. 31125

CAS Registry No.: 2687960-96-7
Formal Name: 4Z,7Z,10Z,13Z,16Z,19Z-docosahexaenoic-21,21,22,22,22-d₅ acid, methyl ester
Synonyms: C22:6 (cis-4,7,10,13,16,19)-d₅ methyl ester, Cervonic Acid-d₅ methyl ester, DHA-d₅ methyl ester, all-cis-4,7,10,13,16,19-DHA-d₅ methyl ester, Methyl all-cis-4,7,10,13,16,19-Docosahexaenoate-d₅, SFE 23:6-d₅
MF: C₂₃H₂₉D₅O₂
FW: 347.6
Chemical Purity: ≥95% (Docosahexaenoic Acid methyl ester)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₅); ≤1% d₀
Supplied as: A 1 mg/ml solution in ethanol
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Docosahexaenoic acid-d₅ methyl ester (DHA-d₅ methyl ester) is intended for use as an internal standard for the quantification of DHA methyl ester (Item No. 10006865) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

DHA-d₅ 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-d₅ methyl ester in these solvents is approximately 100 mg/ml.

Description

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, 06/08/2026

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