

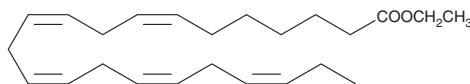
PRODUCT INFORMATION



Docosapentaenoic Acid ethyl ester

Item No. 20669

CAS Registry No.: 119818-40-5
Formal Name: 7Z,10Z,13Z,16Z,19Z-docosapentaenoic acid, ethyl ester
Synonyms: all-*cis*-7,10,13,16,19-Docosapentaenoic Acid ethyl ester, all-*cis*-7,10,13,16,19-DPA ethyl ester, all-*Z*-7,10,13,16,19-DPA ethyl ester, DPA ethyl ester, n-3 DPA ethyl ester, SFE 24:5



MF: C₂₄H₃₈O₂
FW: 358.6
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 ethyl ester is supplied as a solution in ethanol. To change the solvent, simply evaporate the docosapentaenoic acid ethyl ester 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 docosapentaenoic acid ethyl 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 docosapentaenoic acid ethyl ester is needed, it can be prepared by evaporating the ethanol and directly dissolving the neat oil in aqueous buffers. The solubility of docosapentaenoic acid ethyl 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

Docosapentaenoic acid ethyl ester is a derivative of docosapentaenoic acid (DPA; Item No. 90165), an ω-3 fatty acid found in fish oils. Ethyl ester derivatives of fatty acids, including DPA, are often used in formulations designed as dietary supplements.¹⁻³

References

1. James, M.J., Ursin, V.M., and Cleland, L.G. Metabolism of stearidonic acid in human subjects: Comparison with the metabolism of other n-3 fatty acids. *Am. J. Clin. Nutr.* **77**, 1140-1145 (2003).
2. Augustine, A.H., Lowenstein, L.M., Harris, W.S., *et al.* Treatment with ω-3 fatty acid ethyl-ester alters fatty acid composition of lipoproteins in overweight or obese adults with insulin resistance. *Prostaglandins Leukot. Essent. Fatty Acids* **90(2-3)**, 69-75 (2014).
3. Backes, J., Anzalone, D., Hilleman, D., *et al.* The clinical relevance of ω-3 fatty acids in the management of hypertriglyceridemia. *Lipids Health Dis.* **15**, 118-130 (2016).

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