

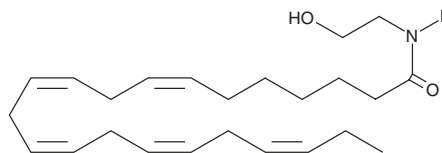
PRODUCT INFORMATION



Docosapentaenoyl Ethanolamide

Item No. 11724

CAS Registry No.: 351317-23-2
Formal Name: N-(2-hydroxyethyl)-7Z,10Z,13Z,16Z,19Z-docosapentaenamide
Synonym: 7(Z),10(Z),13(Z),16(Z),19(Z)-Docosapentaenoyl Ethanolamide
MF: C₂₄H₃₉NO₂
FW: 373.6
Purity: ≥95%
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

Docosapentaenoyl ethanolamide 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 ethanol, DMSO, and dimethyl formamide purged with an inert gas can be used. The solubility of docosapentaenoyl ethanolamide in these solvents is approximately 25, 16, and 20 mg/ml, respectively.

Docosapentaenoyl ethanolamide is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanol solution of docosapentaenoyl ethanolamide should be diluted with the aqueous buffer of choice. Docosapentaenoyl ethanolamide has a solubility of 0.33 mg/ml in a 1:2 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Docosapentaenoyl ethanolamide is a long-chain ω-3 fatty acid with a hydroxyethylamide group attached at the alpha chain.

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, 02/23/2024

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