

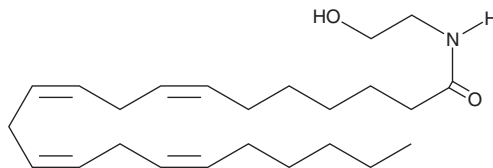
PRODUCT INFORMATION



Docosatetraenoyl Ethanolamide

Item No. 90385

CAS Registry No.: 150314-35-5
Formal Name: N-(2-hydroxyethyl)-7Z,10Z,13Z,16Z-docosatetraenamide
MF: $C_{24}H_{41}NO_2$
FW: 375.6
Purity: $\geq 98\%$
Supplied as: A solution in ethanol
Storage: $-20^{\circ}C$
Stability: ≥ 2 years



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

Laboratory Procedures

Docosatetraenoyl 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 DMSO and dimethyl formamide purged with an inert gas can be used. The solubility of docosatetraenoyl ethanolamide in these solvents is approximately 10 mg/ml.

Docosatetraenoyl ethanolamide is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of docosatetraenoyl ethanolamide should be diluted with the aqueous buffer of choice. Docosatetraenoyl ethanolamide has a solubility of approximately 8 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

Docosatetraenoyl ethanolamide is an endocannabinoid containing docosatetraenoic acid in place of the arachidonate moiety of AEA. Its specific role and relative importance as a cannabinergic neurotransmitter have not been elucidated.¹

Reference

1. Hanus, L., Gopher, A., Almog, S., *et al.* Two new unsaturated fatty acid ethanolamides in brain that bind to the cannabinoid receptor. *J. Med. Chem.* **36**, 3032-3034 (1993).

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, 02/29/2024

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