

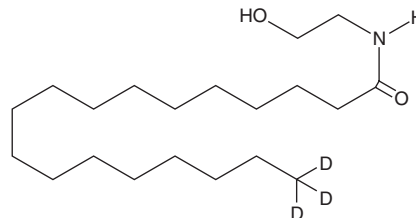
PRODUCT INFORMATION



Stearoyl Ethanolamide-d₃

Item No. 14726

CAS Registry No.: 2749393-37-9
Formal Name: N-(2-hydroxyethyl)-octadecanamide-18,18,18-d₃
Synonyms: Ceramid-d₃, Stearic Acid Ethanolamide-d₃
MF: C₂₀H₃₈D₃NO₂
FW: 330.6
Chemical Purity: ≥98% (Stearoyl Ethanolamide)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₃); ≤1% d₀
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Stearoyl ethanolamide-d₃ is intended for use as an internal standard for the quantification of stearoyl ethanolamide (Item No. 90245) 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).

Stearoyl ethanolamide-d₃ is supplied as a solid. A stock solution may be made by dissolving the stearoyl ethanolamide-d₃ in the solvent of choice, which should be purged with an inert gas. Stearoyl ethanolamide-d₃ is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of stearoyl ethanolamide-d₃ in DMSO is approximately 100 mg/ml and approximately 1 mg/ml in ethanol and DMF.

Description

Stearoyl ethanolamide is a member of the family of fatty N-acyl ethanolamines collectively called anandamides. Stearoyl ethanolamide is the most abundant of several fatty acid ethanolamides produced by the PLD hydrolysis of mouse neuroblastoma cell membrane phospholipids.¹ The specific role of stearoyl ethanolamide in the cannabinergic system is still to be elucidated.

Reference

1. Di Marzo, V., De Petrocellis, L., Sepe, N., *et al.* Biosynthesis of anandamide and related acylethanolamides in mouse J774 macrophages and N18 neuroblastoma cells. *Biochem. J.* **316**(Pt 3), 977-984 (1996).

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, 01/31/2024

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