

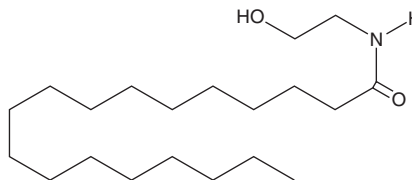
PRODUCT INFORMATION



Stearoyl Ethanolamide

Item No. 90245

CAS Registry No.: 111-57-9
Formal Name: N-(2-hydroxyethyl)-octadecanamide
Synonyms: Ceramid, Stearic Acid Ethanolamide
MF: C₂₀H₄₁NO₂
FW: 327.6
Purity: ≥98%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

Stearoyl ethanolamide is supplied as a crystalline solid. A stock solution may be made by dissolving the stearoyl ethanolamide in the solvent of choice, which should be purged with an inert gas. Stearoyl ethanolamide is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of stearoyl ethanolamide in ethanol and DMF is approximately 2 mg/ml and approximately 100 µg/ml in DMSO.

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. Organic solvent-free aqueous solutions of stearoyl ethanolamide can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of stearoyl ethanolamide in PBS (pH 7.2) is approximately 100 µg/ml. We do not recommend storing the aqueous solution for more than one day.

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, 02/09/2023

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