

PRODUCT INFORMATION



Stearamide

Item No. 21087

CAS Registry No.: 124-26-5
Formal Name: octadecanamide
Synonyms: Amide C18, Amide FA 18:0, NSC 66462, Octadecanamide, Stearic Amide, Stearoyl Amide

MF: $C_{18}H_{37}NO$

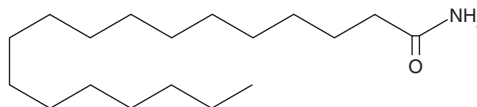
FW: 283.5

Purity: $\geq 98\%$

Supplied as: A crystalline solid

Storage: $-20^{\circ}C$

Stability: ≥ 4 years



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

Laboratory Procedures

Stearamide is supplied as a crystalline solid. A stock solution may be made by dissolving the stearamide in the solvent of choice. Stearamide is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of stearamide in these solvents is approximately 22, 20, and 14 mg/ml, respectively.

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 stearamide can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of stearamide in PBS (pH 7.2) is approximately 50 μ g/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Stearamide is a primary fatty acid amide that is often used in the synthesis of organic chemicals and surfactants. Levels of stearamide are reported to be up-regulated in the serum of patients with hepatic cirrhosis.¹ Thus, it may be a potential biomarker for disordered fatty acid metabolism related to fatty liver diseases.

Reference

1. Lian, J.-S., Liu, W., Hao, S.-R., *et al.* A serum metabonomic study on the difference between alcohol- and HBV-induced liver cirrhosis by ultraperformance liquid chromatography coupled to mass spectrometry plus quadrupole time-of-flight mass spectrometry. *Chin. Med. J. (Engl.)* **124**(9), 1367-1373 (2011).

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, 03/08/2024

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