

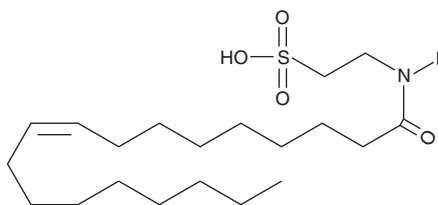
Product Information



N-Oleoyl Taurine

Catalog No. 10005609

CAS Registry No.: 52514-04-2
Formal Name: 2-[(1-oxo-9Z-octadecenyl)amino]-ethanesulfonic acid
MF: C₂₀H₃₉NO₄S
FW: 389.6
Purity: ≥98%
Stability: ≥1 year at -20°C
Supplied as: A solution in DMSO



Laboratory Procedures

For long term storage, we suggest that N-oleoyl taurine be stored as supplied at -20°C. It will be stable for at least one year.

N-Oleoyl taurine is supplied as a solution in DMSO. N-Oleoyl taurine is sparingly soluble in organic solvents such as ethanol and dimethyl formamide.

N-Oleoyl taurine is sparingly soluble in aqueous buffers. Therefore, further dilutions of the organic solvent 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. We do not recommend storing the aqueous solution for more than one day.

Several different arachidonoyl amino acids, including N-arachidonoyl dopamine and N-arachidonoyl serine, have been isolated and characterized from bovine brain.¹ During mass spectral lipidomics analysis of rat brain, a series of fatty acyl amides of a third amino acid, taurine, were discovered.² This novel class of compounds is present in kidney and activates members of the transient receptor potential (TRP) family of calcium channels.³ N-Oleoyl taurine is an amino-acyl endocannabinoid isolated from rat brain that may activate TRPV1 and TRPV4.³

References

- Huang, S.M., Bisogno, T., Petros, T.J., *et al.* Identification of a new class of molecules, the arachidonoyl amino acids, and characterization of one member that inhibits pain. *J. Biol. Chem.* **276(46)**, 42639-42644 (2001).
- Cravatt, B.F. and Maxey, K.M. Personal communication. (2004).
- Saghatelian, A., McKinney, M.K., Bandell, M., *et al.* A FAAH-regulated class of N-acyl taurines that activates TRP ion channels. *Biochemistry* **45(30)**, 9007-9015 (2006).

Related Products

1-Arachidonoyl Glycerol - Cat. No. 62150 • 2-Arachidonoyl Glycerol - Cat. No. 62160 • Arachidonoyl Ethanolamide - Cat. No. 90050 • Arachidonoyl 2-Fluoroethylamide - Cat. No. 90054 • O-Arachidonoyl Ethanolamine (hydrochloride) - Cat. No. 91050 • N-Arachidonoyl Taurine - Cat. No. 10005537 • N-Stearoyl Taurine - Cat. No. 10005610 • N-Palmitoyl Taurine - Cat. No. 10005611

WARNING: THIS PRODUCT IS NOT FOR HUMAN OR ANIMAL DISEASE DIAGNOSIS OR THERAPEUTIC DRUG USE.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent under separate cover to the MSD supervisor at your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will meet our specifications at the time of delivery.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a refund of the purchase price, or at Cayman's option, the replacement, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy** located on our website and in our catalog.

Copyright Cayman Chemical Company, 08/01/2008

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com