

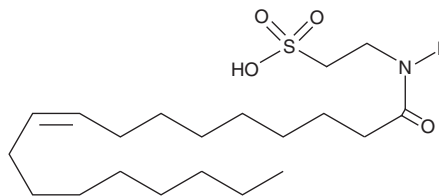
PRODUCT INFORMATION



N-Oleoyl Taurine

Item 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%
Supplied as: A solution in DMSO
Storage: -20°C
Stability: ≥1 year



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

Description

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

1. 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).
2. Cravatt, B.F. and Maxey, K.M. Personal communication. (2004).
3. 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).

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, 06/03/2020

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