

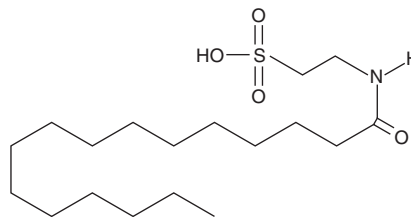
PRODUCT INFORMATION



N-Palmitoyl Taurine

Item No. 10005611

CAS Registry No.: 83982-06-3
Formal Name: 2-[(1-oxohexadecyl)amino]-ethanesulfonic acid
MF: C₁₈H₃₇NO₄S
FW: 363.6
Purity: ≥98%
UV/Vis.: λ_{max}: 203, 256 nm
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.

Description

Several different arachidonoyl amino acids, including N-arachidonoyl dopamine (NADA) and N-arachidonoyl serine (ARA-S), have been isolated and characterized from bovine brain.¹ During mass spectral lipidomic analysis of rat brain, a series of fatty acyl amides of a third amino acid, taurine, was discovered. N-Palmitoyl taurine is a prominent amino-acyl endocannabinoid isolated from rat brain during lipidomics profiling. Its function is currently under investigation.

Reference

1. Greig, G.M., Francis, D.A., Falgoutyret, J.P., *et al.* The interaction of arginine 106 of human prostaglandin G/H synthase-2 with inhibitors is not a universal component of inhibition mediated by nonsteroidal anti-inflammatory drugs. *Mol. Pharmacol.* **52(5)**, 829-838 (1997).

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, 11/10/2023

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