

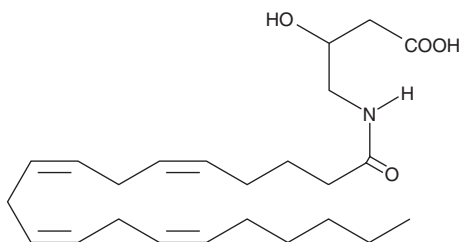
PRODUCT INFORMATION



N-Arachidonoyl-3-hydroxy-γ-Aminobutyric Acid

Item No. 10158

CAS Registry No.: 959761-62-7
Formal Name: 4-[[[(3-hydroxy)-1-oxo-5Z,8Z,11Z,14Z-eicosatetraenyl]amino]-butanoic acid
Synonym: NAG-3H-ABA
MF: C₂₄H₃₉NO₄
FW: 405.6
Purity: ≥98%
Stability: ≥1 year at -20°C
Supplied as: A solution in methyl acetate



Laboratory Procedures

For long term storage, we suggest that N-arachidonoyl-3-hydroxy-γ-aminobutyric acid (NAG-3H-ABA) be stored as supplied at -20°C. It should be stable for at least one year.

NAG-3H-ABA is supplied as a solution in methyl acetate. To change the solvent, simply evaporate the methyl acetate under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide purged with an inert gas can be used. The solubility of NAG-3H-ABA in these solvents is approximately 30, 15, and 20 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. If an organic solvent-free solution of NAG-3H-ABA is needed, it can be prepared by evaporating the ethanol and directly dissolving the neat oil in aqueous buffers. The solubility of NAG-3H-ABA in PBS (pH 7.2) is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Several different arachidonoyl amino acids, including NAG-3H-ABA, have been isolated and characterized from bovine brain.¹ The glycine congener (NAGly; Item No. 90051) was further characterized and found to suppress formalin-induced pain in rats. NAG-3H-ABA was also found in rat brain by LC-MS techniques, but has not been fully characterized to date. Most arachidonoyl amino acids are poor ligands for the CB₁ receptor.

Reference

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).

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, 08/13/2015

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