

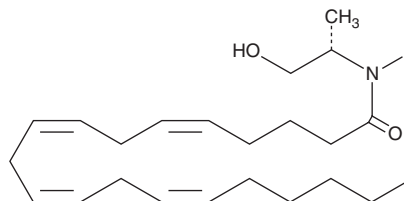
PRODUCT INFORMATION



S-1 Methanandamide

Item No. 90072

CAS Registry No.: 157182-50-8
Formal Name: N-(2-hydroxy-1S-methylethyl)-5Z,8Z,11Z,14Z-eicosatetraenamide
Synonym: (S)-(-)-Arachidonyl-1'-Hydroxy-2'-Propylamide
MF: C₂₃H₃₉NO₂
FW: 361.6
Purity: ≥98%
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥2 years



Special Conditions: Oxygen and light sensitive

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

Laboratory Procedures

S-1 Methanandamide is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as DMSO, dimethyl formamide, or acetonitrile purged with an inert gas or nitrogen can be used.

S-1 Methanandamide is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of S-1 methanandamide should be diluted with the aqueous buffer of choice. We do not recommend storing the aqueous solution for more than one day.

Description

S-1 Methanandamide is a CB₁ receptor ligand but is less potent than the C-1 (R) methyl isomer. It inhibits electrically evoked contractions in isolated mouse vasa deferentia with an IC₅₀ of 230 nM. The binding affinity of S-1 methanandamide for CB₁ receptors is less than that of AEA, with a K_i of 175 nM for the displacement of radiolabeled CP 55,940.¹

Reference

1. Abadji, V., Lin, S., Taha, G., *et al.* (R)-Methanandamide: A chiral novel anandamide possessing higher potency and metabolic stability. *J. Med. Chem.* **37**, 1889-1893 (1994).

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, 02/28/2024

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