

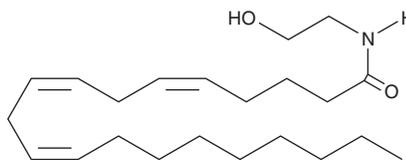
PRODUCT INFORMATION



5(Z),8(Z),11(Z)-Eicosatrienoic Acid Ethanolamide

Item No. 90195

CAS Registry No.: 169232-04-6
Formal Name: N-(2-hydroxyethyl)-5Z,8Z,11Z-eicosatrienamide
Synonym: Mead Acid Ethanolamide
MF: C₂₂H₃₉NO₂
FW: 349.6
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

Mead acid ethanolamide 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 and dimethyl formamide purged with an inert gas can be used. The solubility of mead acid ethanolamide in these solvents is approximately 10 and 30 mg/ml, respectively.

Mead acid ethanolamide is a lipid soluble compound, and therefore is not directly soluble in aqueous solutions. 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.

Description

Mead acid ethanolamide is essentially identical to AEA in its agonist binding to CB₁ and CB₂ receptors. In L cells expressing the human CB₁ receptor, the K_i value for mead acid ethanolamide and AEA binding is 753 nM. In AtT-20 cells expressing the human CB₂ receptor, mead acid ethanolamide and AEA bind with a K_i value of 1,810 nM.¹

Reference

1. Priller, J., Briley, E.M., Mansouri, J., *et al.* Mead ethanolamide, a novel eicosanoid, is an agonist for the central (CB₁) and peripheral (CB₂) cannabinoid receptors. *Mol. Pharmacol.* **48**, 288-292 (1995).

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, 02/29/2024

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