

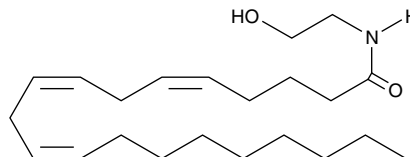
Product Information



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%
Stability: ≥1 year at -20°C
Supplied as: A solution in ethanol
Misc.: Oxygen and light sensitive



Laboratory Procedures

For long term storage, we suggest that mead acid ethanolamide be stored as supplied at -20°C. It should be stable for at least one year.

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.

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 ArT-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).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/90195

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY. NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. 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

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery**.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees. Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our Warranty and Limitation of Remedy located on our website and in our catalog.

Copyright Cayman Chemical Company, 02/05/2014

Cayman Chemical

Mailing address
1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone
(800) 364-9897
(734) 971-3335

Fax
(734) 971-3640

E-Mail
custserv@caymanchem.com

Web
www.caymanchem.com