

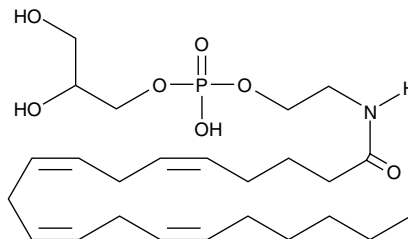
Product Information



Glycerophospho-N-Arachidonoyl Ethanolamine

Item No. 10011347

CAS Registry No.: 201738-25-2
Formal Name: mono(2,3-dihydroxypropyl)-
mono[2-[[[(5Z,8Z,11Z,14Z)-1-oxo-
5,8,11,14-eicosatetraen-1-yl]ethyl]ester
phosphoric acid
Synonyms: Glycerophosphoanandamide,
Glycerophospho-Arachidonoyl
Ethanolamide, GP-NArE
MF: C₂₅H₄₄NO₇P
FW: 501.6
Purity: ≥97%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid



Laboratory Procedures

For long term storage, we suggest that glycerophospho-N-arachidonoyl ethanolamine (GP-NArE) be stored as supplied at -20°C. It should be stable for at least two years.

GP-NArE is supplied as a crystalline solid. A stock solution may be made by dissolving the GP-NArE in an organic solvent purged with an inert gas. GP-NArE is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of GP-NArE in these solvents is approximately 20 mg/ml.

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. Organic solvent-free aqueous solutions of GP-NArE can be prepared by directly dissolving the crystalline compound in aqueous buffers. The solubility of GP-NArE in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

N-Acylated ethanolamines (NAE) are naturally-occurring lipids that have diverse bioactivities. The different types of NAE can be derived from glycerophospho-linked precursors by the activity of glycerophosphodiesterase 1 (GDE₁).¹ GP-NArE is the precursor of arachidonoyl ethanolamide (AEA), also known as anandamide. AEA is an endogenous cannabinoid neurotransmitter that binds to both central cannabinoid (CB₁) and peripheral cannabinoid (CB₂) receptors.² It inhibits the specific binding of [³H]-HU-243 to synaptosomal membranes with a K_i value of 52 nM, compared to 46 nM for Δ⁹-THC.³

References

1. Simon, G.M. and Cravatt, B.F. Anandamide biosynthesis catalyzed by the phosphodiesterase GDE₁ and detection of glycerophospho-N-acyl ethanolamine precursors in mouse brain, [In Press] *J. Biol. Chem.* (2008).
2. Felder, C.C., Briley, E.M., Axelrod, J., *et al.* Anandamide, an endogenous cannabimimetic eicosanoid, binds to the cloned human cannabinoid receptor and stimulates receptor-mediated signal transduction. *Proc. Natl. Acad. Sci. USA* **90**, 7656-7660 (1993).
3. Devane, W.A., Hanus, L., Breuer, A., *et al.* Isolation and structure of a brain constituent that binds to the cannabinoid receptor. *Science* **258**, 1946-1949 (1992).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/10011347

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, 06/24/2014

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