

PRODUCT INFORMATION

Glycerophospho-N-Eicosapentaenoyl Ethanolamine

Item No. 90001842

Formal Name: 2,3-dihydroxypropyl
(2-((5Z,8Z,11Z,14Z,17Z)-icosa-
5,8,11,14,17-pentaenamido)ethyl)
hydrogen phosphate

Synonym: GP-NEPEA

MF: C₂₅H₄₂NO₇P

FW: 499.6

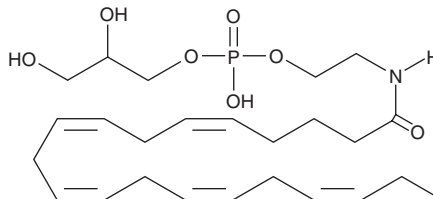
Purity: ≥95%

UV/Vis.: λ_{max}: 209 nm

Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥4 years



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

Laboratory Procedures

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

GP-NEPEA is supplied as a crystalline solid. A stock solution may be made by dissolving the GP-NEPEA in the solvent of choice. GP-NEPEA is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of GP-NEPEA 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-NEPEA can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of GP-NEPEA in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

N-Acylated ethanolamines (NAEs) are naturally-occurring lipids that have diverse bioactivities. For example, arachidonoyl ethanolamide (Item No. 90050) is an endogenous neurotransmitter that evokes cellular responses by activating the cannabinoid (CB) receptors, central CB₁ and peripheral CB₂. The different types of NAEs are derived from glycerophospho-linked precursors by the activity of glycerophosphodiesterase 1.¹ GP-NEPEA is the metabolic precursor of eicosapentaenoyl ethanolamide (EPEA; Item No. 10964). EPEA is an NAE that inhibits dietary-restriction-induced lifespan extension in wild type and TOR pathway mutant nematodes.² EPEA also has anti-inflammatory properties, suppressing the expression of IL-6 and MCP-1 in 3T3-L1 adipocytes in response to lipopolysaccharide.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. *J. Biol. Chem.* **283**, 9341-9349 (2008).
2. Lucanic, M., Held, J.M., Vantipalli, M.C., et al. N-acyl ethanolamine signalling mediates the effect of diet on lifespan in *Caenorhabditis elegans*. *Nature* **473**(7346), 226-9 (2011).
3. Balvers, M.G., Verhoeckx, K.C., Plastina, P., et al. Docosahexaenoic acid and eicosapentaenoic acid are

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, 04/09/2024

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