

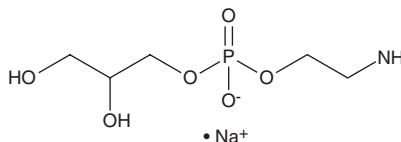
PRODUCT INFORMATION



Glycerophosphorylethanolamine (sodium salt)

Item No. 34460

CAS Registry No.: 883288-78-6
Formal Name: phosphoric acid, mono(2-aminoethyl) mono(2,3-dihydroxypropyl) ester, monosodium salt
Synonyms: Glycerol 3-phosphorylethanolamine, Glycerolphosphorylethanolamine
MF: C₅H₁₃NO₆P • Na
FW: 237.1
Purity: ≥90%
Supplied as: A 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

Glycerophosphorylethanolamine (sodium salt) is supplied as a solid. A stock solution may be made by dissolving the glycerophosphorylethanolamine (sodium salt) in water. Glycerophosphorylethanolamine (sodium salt) is slightly soluble in water. We do not recommend storing the aqueous solution for more than one day.

Description

Glycerophosphorylethanolamine is an active phosphodiester metabolite of phosphatidylethanolamine.^{1,2} It promotes aggregation of amyloid-β (1-40) (Aβ40) *in vitro*, and levels of glycerophosphorylethanolamine are elevated in postmortem brains isolated from patients with Alzheimer's disease.

References

1. Klunk, W.E., Xu, C.J., McClure, R.J., *et al.* Aggregation of β-amyloid peptide is promoted by membrane phospholipid metabolites elevated in Alzheimer's disease brain. *J. Neurochem.* **69(1)**, 266-272 (1997).
2. Blusztajn, J.K., Lopez Gonzalez-Coviella, I., Logue, M., *et al.* Levels of phospholipid catabolic intermediates, glycerophosphocholine and glycerophosphoethanolamine, are elevated in brains of Alzheimer's disease but not of Down's syndrome patients. *Brain Res.* **536(1-2)**, 240-244 (1990).

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, 11/29/2022

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