

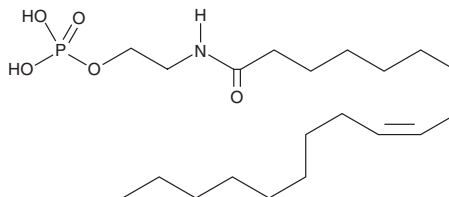
PRODUCT INFORMATION



Oleoyl Ethanolamide Phosphate

Item No. 22886

CAS Registry No.: 24435-25-4
Formal Name: N-[2-(phosphonooxy)ethyl]-9Z-octadecenamide
Synonyms: NAEPA, N-oleoyl ethanolamide phosphoric acid, OEA-P
MF: C₂₀H₄₀NO₅P
FW: 405.5
Purity: ≥95%
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

Oleoyl ethanolamide phosphate (OEA-P) is supplied as a crystalline solid. A stock solution may be made by dissolving the OEA-P in the solvent of choice, which should be purged with an inert gas. OEA-P is soluble in DMSO.

Description

OEA-P is a lysophosphatidic acid (LPA) mimetic, LPA receptor agonist, and phosphate ester of oleoyl ethanolamide (OEA; Item No. 90265).^{1,2} It selectively increases [³⁵S]GTPγS binding to HEK293T cell membranes expressing LPA₁ and LPA₂ over LPA₃ receptors.² OEA-P induces calcium mobilization and inhibits forskolin-induced cAMP accumulation in MDA-MB-231 cells (EC₅₀s = 1.2 and 101 nM, respectively).¹

References

1. Lynch, K.R., Hopper, D.W., Carlisle, S.J., *et al.* Structure/activity relationships in lysophosphatidic acid: The 2-hydroxyl moiety. *Mol. Pharmacol.* **52**(1), 75-81 (1997).
2. Im, D.S., Heise, C.E., Harding, M.A., *et al.* Molecular cloning and characterization of a lysophosphatidic acid receptor, Edg-7, expressed in prostate. *Mol. Pharmacol.* **57**(4), 753-759 (2000).

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, 12/01/2022

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