PRODUCT INFORMATION



1-Stearoyl-2-Adrenoyl-sn-glycero-3-PE

Item No. 33661

| CAS Registry No.: Formal Name: | 131350-53-3 (7Z,10Z,13Z,16Z)-7,10,13,16-docosatetraenoic acid (1R)-1-[[[(2-aminoethoxy)hydroxyphosphinyl]oxy] methyl]-2-[(1-oxooctadecyl)oxy]ethyl ester | |
|-----------------------------------|--|--|
| Synonyms: | 1-Octadecanoyl-2-Docosatetraenoyl- sn-glycero-3-Phosphatidylethanolamine, 1-Octadecanoyl-2-Docosatetraenoyl-sn-glycero- 3-Phosphoethanolamine, 1-Octadecanoyl-2- Docosatetraenoyl-sn-glycero-3-PE, 1-Stearoyl-2- Docosatetranenoyl-sn-glycero-3-PE, PE(18:0/22:4) | |
| MF: | C ₄₅ H ₈₂ NO ₈ P | ОН |
| FW: | 796.1 | |
| Purity: | ≥95% | |
| Supplied as: | A solution in chloroform | |
| Storage: | -20°C | |
| Stability: | ≥1 year | |
| Information represents | the product specifications. Batch specific analytical results are | provided on each certificate of analysis |

Description

1-Stearoyl-2-adrenoyl-sn-glycero-3-PE is a phospholipid that contains stearic acid (Item No. 10011298) and adrenic acid (Item No. 90300) at the sn-1 and sn-2 positions, respectively. 1-Stearoyl-2-adrenoyl-snglycero-3-PE levels are inversely correlated with subject age in mitochondria isolated from human postmortem hippocampus.¹

Reference

1. Hancock, S.E., Friedrich, M.G., Mitchell, T.W., et al. Decreases in phospholipids containing adrenic and arachidonic acids occur in the human hippocampus over the adult lifespan. Lipids 50(9), 861-872 (2015).

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.

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, 05/05/2021

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