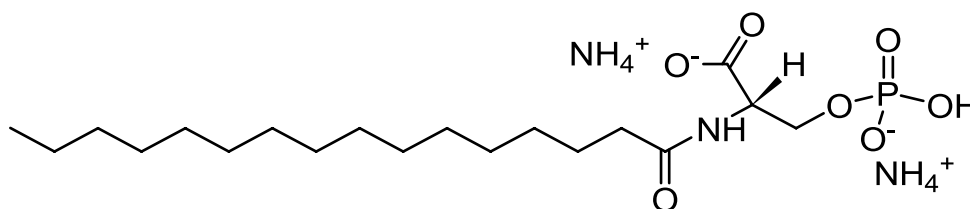


TECHNICAL DATA SHEET

N-Palmitoyl-Serine-Phosphoric Acid

Catalog Number	800715	Physical state	Powder
Purity	> 99%	Transition temp.	No data
CAS	799268-44-3	CMC	No data
Synonyms	N-16:0-Serine-PO ₄	PKA	No data
Molec. Formula	C ₁₉ H ₄₄ N ₃ O ₇ P	TLC mobile phase	C:M:W & C:M:A *, 65:35:8, v/v
MW	457.550	Exact Mass	457.292
Percent composition	C 49.88% H 9.69% N 9.18% O 24.48% P 6.77%		
Stability	Store in <-20°C freezer for up to one year		
Solubility	See note in Product Use section below		
Web link	800715		

*chloroform:methanol:water and chloroform:methanol:acetone



Description:

Lysophosphatidic acid (LPA) receptor modulators include N-palmitoyl serine phosphoric acid and N-palmitoyl-tyrosine phosphoric acid. N-palmitoyl serine phosphoric acid is a competitive inhibitor of the LPA receptor in *Xenopus* oocytes (Neidlinger et al, 2006; Liliom et al, 1996). However, in mammalian cells, N-palmitoyl-serine phosphoric acid may act as an agonist for the LPA receptor (Hooks et al, 1998; An et al, 1998). LPA is a lipid mediator that acts similar to growth factors through G-protein coupled plasma membrane receptors (Neidlinger et al, 2006; Liliom et al, 1996; Hooks et al, 1998; An et al, 1998; Jan et al, 2003). LPA may play a role in platelet aggregation, smooth muscle contraction, vasoactive changes, cytoskeletal reorganization and cell proliferation (Neidlinger et al, 2006). N-palmitoyl serine phosphoric acid may also play a role in signal transduction in mammalian cells by increasing intracellular calcium (An et al, 1998; Jan et al, 2003).

Product use:

A range of 0.1-20 µM of N-palmitoyl serine phosphoric acid can be used for cell studies. Make a stock solution of 5 mM in DMSO and store at -20°C. Diluted N-palmitoyl serine phosphoric acid can be directly added to the study medium (Jan et al, 2003). An et al. dissolved N-palmitoyl-serine and N-palmitoyl-tyrosine phosphoric acid in 0.1 mL PBS containing 0.1 mg/mL human serum albumin before adding to cells (An et al, 1998). In *X.laevis* studies, these LPA inhibitors were dissolved in DMSO at 1 mM and filtered through a 0.45 µM membrane filter before injection (Liliom et al, 1996).

References:

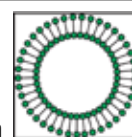
- Khandoga et al (2008) Lysophosphatidic acid-induced platelet shape change revealed through LPA(1-5)receptor-selective probes and albumin. Platelets 19(6):415-27
- Neidlinger NA et al (2006) Hydrolysis of phosphatidylserine-exposing red blood cells by secretory phospholipase A2 generates lysophosphatidic acid and results in vascular dysfunction. J Biol Chem. 281(2): 775-81
- Jan CR et al (2003) Novel effect of N-palmitoyl-L-serine phosphoric acid on cytosolic Ca²⁺ levels in human osteoblasts. Pharmacol Toxicol 93(2): 71-6
- Hooks SB et al (1998) Characterization of a receptor subtype-selective lysophosphatidic mimetic. Molec Pharm. 53: 188-94
- An S et al (1998) Recombinate human G protein-coupled lysophosphatidic receptors mediate intracellular calcium mobilization. Molec Pharm. 54: 881-8
- Liliom K et al (1996) N-Palmitoyl-serine and N-palmitoyl-tyrosine phosphoric acids are selective competitive antagonists of the lysophosphatidic acid receptors. Molec Pharm 50:616-23

Related products: [N-palmitoyl-tyrosine phosphoric acid](#)
[Receptor agonist/antagonist](#)

MSDS: Available on Avanti's website for product number 800715

Avanti Polar Lipids, Inc., 700 Industrial Park Drive, Alabaster, AL 35007-9105, U.S.A.

- (800) 227-0651 • (205) 663-2494 • Fax (800) 229-1004 • Fax (205) 663-0756
- Email Orders: orders@avantilipids.com • Email Inquiries: info@avantilipids.com
- Email Technical Questions: technical@avantilipids.com • Visit www.avantilipids.com



Avanti[®]
POLAR LIPIDS, INC.