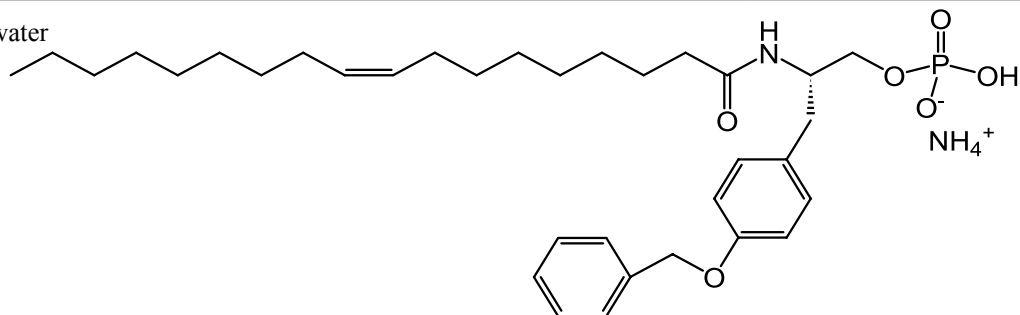


TECHNICAL DATA SHEET

(S)-Phosphoric acid mono-[3-(4-benzyloxy-phenyl)-2-octadec-9-enoylamino-propyl] ester (ammonium salt) (VPC 12249(S))

Catalog Number	857341	Physical state	Powder
Purity	> 99%	Transition temp.	No data
CAS	799268-73-8	CMC	No data
Synonyms	LPA ₁ /LPA ₃ receptor antagonist; VPC 12249	pK_a	No data
Molec. Formula	C ₃₄ H ₅₅ N ₂ O ₆ P	TLC mobile phase	C:M:W*, 65:35:8, v/v
MW	618.784	Exact Mass	618.380
Percent composition	C 65.99% H 8.96% N 4.53% O 15.51% P 5.01%		
Stability	Store in <-20°C freezer for up to 6 months. Aliquot suspensions (1 mM) and store frozen.		
Solubility	Suspend VPC 12249 in 3% BSA (fatty acid free Bovine Serum Albumin) in water at a lipid concentration of 1 mM.		
Web link	857341		

*chloroform:methanol:water



Description:

Lysophospholipids play a role in a broad spectrum of cellular functions, including signal transduction, membrane trafficking and cell growth, migration and survival (Sigal *et al*, 2005). The actions of lysophospholipids, including lysophosphatidic acid (LPA) and sphingosine 1-phosphate (S1P), have been studied through specific interactions with ten G-protein-coupled receptors (LPA₁₋₅ and S1P₁₋₅) (Skoura and Hla, 2009) and with the nuclear receptor PPAR- γ (peroxisome-proliferator-activated receptor- γ) (Prestwich *et al*, 2005). By defining specific receptor agonists and antagonists, lysophospholipids have been implicated in such diverse pathophysiological states such as cancer, autoimmune diseases, atherosclerosis (Gardell *et al*, 2006; Prestwich *et al*, 2005), immunodeficiency, ischemia-reperfusion injury (Prestwich *et al*, 2005), neuropathic pain and obesity (Gardell *et al*, 2006). Therefore lysophospholipid receptors have emerged as drug targets for therapeutic intervention (Gardell *et al*, 2006).

VPC 12249(S) is a LPA₁ and LPA₃ receptor antagonist.

How to use:

Please use the following web links for [TLC](#) or [liposome preparation](#)

References:

- Skoura A, Hla T (2009) Lysophospholipid receptors in vertebrate development, physiology, and pathology. *J Lipid Res.* 2009 Apr;50 Suppl:S293-8
- Gardell SE, Dubin AE, Chun J (2006) Emerging medicinal roles for lysophospholipid signaling. *Trends Molec Med* 12(2): 65-75
- Sigal YJ, McDermott MI, Morris AJ (2005) Integral membrane lipid phosphatases/phosphotransferases: common structure and diverse functions. *Biochem J* 387: 281-293
- Chun, J (2005) Lysophospholipids in the nervous system. *Prostaglandins & other Lipid Mediators* 77: 46-51
- Prestwich GD *et al* (2005) New metabolically stabilized analogues of lysophosphatidic acid: agonists, antagonists and enzyme inhibitors. *Biochem Soc Trans.* 33: 1357-1361
- Davis MD *et al* (2005) Spingosine-1-phosphate analogs as receptor antagonists. *J Biol Chem* 280(11): 9833-9841
- Santos WL *et al* (2004) Synthesis and biological evaluation of phosphonic and thiophosphoric acid derivatives of lysophosphatidic acid. *Bioorg Med Chem Lett* 14:3473-3476.

Related products: [Receptor Agonist/Antagonist](#)

MSDS: Available at www.avantilipids.com for Product Number 857341

Avanti Polar Lipids, Inc., 700 Industrial Park Drive, Alabaster, AL 35007-9105, U.S.A.
Phone • (800) 227-0651 • (205) 663-2494 • Fax (800) 229-1004 • (205) 663-0756
Email • orders@avantilipids.com • Technical Questions: technical@avantilipids.com
• Inquiries: info@avantilipids.com • Analytical: analytical@avantilipids.com

