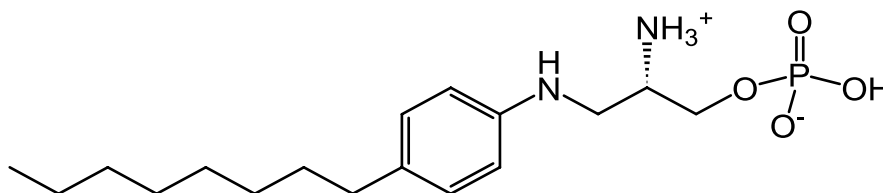


# TECHNICAL DATA SHEET

## (S)-Phosphoric acid mono-[2-amino-3-(4-octyl-phenylamino)-propyl] ester (VPC 24191)

<b>Catalog Number</b>	857365	<b>Physical state</b>	Powder
<b>Purity</b>	> 99%	<b>Transition temp.</b>	No data
<b>CAS</b>	799268-83-0	<b>CMC</b>	No data
<b>Synonyms</b>	S1P <sub>1</sub> /S1P <sub>3</sub> competitive agonist; VPC 24191	<b>pK<sub>a</sub></b>	No data
<b>Molec. Formula</b>	C <sub>17</sub> H <sub>31</sub> N <sub>2</sub> O <sub>4</sub> P	<b>TLC mobile phase</b>	C:M:W*, 65:35:8, v/v
<b>MW</b>	358.413	<b>Exact Mass</b>	358.202
<b>Percent composition</b>	C 56.97% H 8.72% N 7.82% O 17.86% P 8.64%		
<b>Stability</b>	Store in <-20°C freezer for up to 6 months		
<b>Solubility</b>	Dissolve to 20mM in DMSO/1N HCl (95:5 v/v). Dilute (1:20) immediately into 3% aqueous fatty acid free BSA. Final stock is 1mM lipid, 95 parts BSA, 5 parts acidified DMSO. Aliquot and store at <-20°C; avoid freeze/thaw		
<b>Web link</b>	<a href="#">857365</a>		

\*chloroform:methanol:water



### Description:

Sphingosine-1-phosphate (S1P) is a lysophospholipid mediator that evokes a variety of cellular responses by stimulation of five members of the endothelial cell differentiation gene receptor family. The endothelial cell differentiation gene receptors are G-protein coupled receptors that, upon stimulation, propagate second messenger signals via activation of heterotrimeric G-protein subunits and dimers. Ultimately, this S1P-driven signaling results in cell survival, increased cell migration, and, often, mitogenesis. (Davis *et al*, 2005).

VPC 24191 is a competitive agonist at the S1P<sub>1</sub> and S1P<sub>3</sub> receptors.

### 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
- Davis MD *et al* (2005) Sphingosine-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](http://www.avantilipids.com) for Product Number 857365