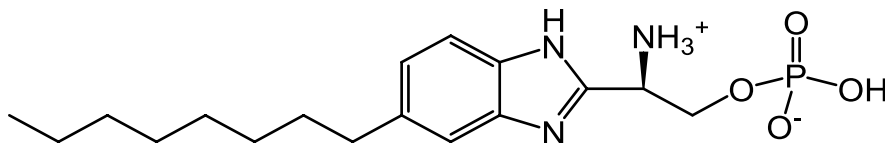


# TECHNICAL DATA SHEET

## (R)-Phosphoric acid mono-[2-amino-2-(6-octyl-1H-benzimidazol-2-yl) ethyl] ester (VPC 23153)

Catalog Number	857367	Physical state	Powder
Purity	> 99%	Transition temp.	No data
CAS	787582-98-3	CMC	No data
Synonyms	S1P <sub>4</sub> competitive agonist; VPC 23153	pK <sub>a</sub>	No data
Molec. Formula	C <sub>17</sub> H <sub>28</sub> N <sub>3</sub> O <sub>4</sub> P	TLC mobile phase	C:M:W*, 65:35:8, v/v
MW	369.396	Exact Mass	369.182
Percent composition	C 52.27% H 7.64% N 11.38% O 17.32% P 8.38%		
Stability	Store in <-20°C freezer for up to 6 months		
Solubility	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		
Web link	<a href="#">857367</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 23153 is a competitive agonist at the S1P<sub>4</sub> receptor.

### 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) Spingosine-1-phosphate analogs as receptor antagonists. *J Biol Chem* 280(11): 9833-9841
- Clemens, J.J., Davis, M.D., Lynch, K.R., Macdonald, T.L. (2004) Synthesis of benzimidazole based analogues of sphingosine-1-phosphate: discovery of potent, subtype-selective S1P<sub>4</sub> receptor agonists. *Bioorg Med Chem Lett* 14:4903-6.

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

**MSDS:** Available at [www.avantilipids.com](http://www.avantilipids.com) for Product Number 857367