

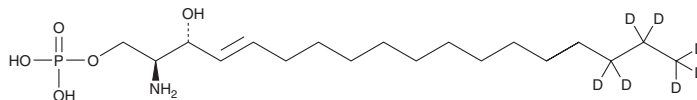
# PRODUCT INFORMATION



## Sphingosine-1-phosphate-d<sub>7</sub> (d18:1)

Item No. 10008121

**CAS Registry No.:** 2260670-15-1  
**Formal Name:** (2S,3R,4E)-2-amino-4-octadecene-16,16,17,17,18,18,18-d<sub>7</sub>-1,3-diol, 1-(dihydrogen phosphate)  
**Synonyms:** S1P-d<sub>7</sub> (d18:1), Sphingosine-1-Phosphoric Acid-d<sub>7</sub>  
**MF:** C<sub>18</sub>H<sub>31</sub>D<sub>7</sub>NO<sub>5</sub>P  
**FW:** 386.5  
**Chemical Purity:** ≥95% (Sphingosine-1-phosphate (d18:1))  
**Deuterium Incorporation:** ≥99% deuterated forms (d<sub>1</sub>-d<sub>7</sub>); ≤1% d<sub>0</sub>  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

Sphingosine-1-phosphate-d<sub>7</sub> (d18:1) is intended for use as an internal standard for the quantification of sphingosine-1-phosphate (S1P; Item No. 62570) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

### Description

S1P is the product of phosphorylation of sphingosine by sphingosine kinase that is secreted from cells and acts as an agonist at S1P receptors.<sup>1</sup> It increases intracellular calcium levels in TAg-Jurkat cells expressing S1P<sub>1</sub> and G<sub>q/15</sub>, which allows for phospholipase C stimulation by G<sub>i</sub> proteins, when used at a concentration of 200 nM, as well as in TAg-Jurkat cells expressing S1P<sub>2</sub> and S1P<sub>3</sub> receptors (EC<sub>50</sub>s = 8 and 11 nM, respectively).<sup>2</sup> Intra- and extracellular levels of S1P vary spatially allowing it to function as an autocrine or paracrine factor, respectively, and dysregulation of S1P levels are associated with various disease states, such as inflammation and autoimmunity.<sup>3</sup> S1P has a wide variety of effects, including an involvement in cell growth, angiogenesis, immunity, and neuroprotection.

### References

1. Sanchez, T. and Hla, T. Structural and functional characteristics of S1P receptors. *J. Cell. Biochem.* **92**(5), 913-922 (2004).
2. An, S., Bleu, T., and Zheng, Y. Transduction of intracellular calcium signals through G protein-mediated activation of phospholipase C by recombinant sphingosine 1-phosphate receptors. *Mol. Pharmacol.* **55**(5), 787-794 (1999).
3. Cartier, A. and Hla, T. Sphingosine 1-phosphate: Lipid signaling in pathology and therapy. *Science* **366**(6463), (2019).

#### 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.

#### WARRANTY AND LIMITATION OF REMEDY

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, 01/22/2024

#### 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](http://WWW.CAYMANCHEM.COM)