

PRODUCT INFORMATION



Sphingosine-1-phosphate (d17:1)

Item No. 22498

CAS Registry No.: 474923-27-8

Formal Name: 2S-amino-4E-heptadecene-1,3R-diol,
1-(dihydrogen phosphate)

Synonyms: D-erythro-Sphingosine C17-phosphate,
S1P (d17:1)

MF: $C_{17}H_{36}NO_5P$

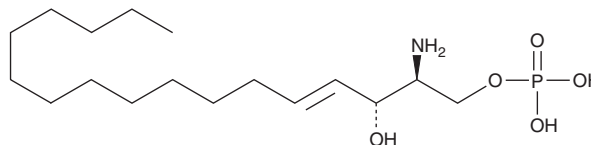
FW: 365.5

Purity: $\geq 95\%$

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 (S1P) (d17:1) is supplied as a crystalline solid. A stock solution may be made by dissolving the S1P (d17:1) in the solvent of choice. S1P (d17:1) is soluble in 0.3 M sodium hydroxide, which should be purged with an inert gas, at a concentration of approximately 4 mg/ml.

Description

S1P (d17:1) is a synthetic sphingolipid that has been used as an internal standard for the quantification of S1P (d18:1) (Item No. 62570) in plasma and tissue samples.^{1,2}

References

1. Qu, F., Wu, C.S., Hou, J.F., *et al.* Sphingolipids as new biomarkers for assessment of delayed-type hypersensitivity and response to triptolide. *PLoS One* **7**(12), e52454 (2012).
2. Berdyshev, E.V., Gorshkova, I.A., Garcia, J.G., *et al.* Quantitative analysis of sphingoid base-1-phosphates as bisacetylated derivatives by liquid chromatography-tandem mass spectrometry. *Anal. Biochem.* **339**(1), 129-136 (2005).

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, 02/15/2024

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