

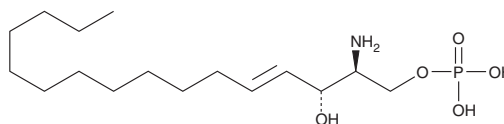
PRODUCT INFORMATION



Sphingosine-1-phosphate (d16:1)

Item No. 9002921

CAS Registry No.: 709026-60-8
Formal Name: 2S-amino-1-(dihydrogen phosphate),
4E-hexadecene-1,3R-diol
Synonyms: C16 S1P, C16 Sphingosine-1-
phosphate, S1P (d16:1)
MF: C₁₆H₃₄NO₅P
FW: 351.4
Purity: ≥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

C16 Sphingosine-1-phosphate (C16 S1P) is supplied as a crystalline solid. C16 S1P is sparingly soluble in organic solvents. Aqueous solutions can be prepared by directly dissolving the crystalline solid in basic buffers. The solubility of C16 S1P in 0.3 M NaOH is approximately 4 mg/ml.

Description

C16 S1P is a derivative of sphingosine-1-phosphate (S1P; Item No. 62570) that binds to S1P₁/EDG-1, S1P₃/EDG-3, and S1P₂/EDG-5 receptors with affinities of 115%, 83%, and 103%, respectively, relative to S1P in CHO cells.¹ C16 S1P was increased in postmortem primary open angle glaucoma trabecular meshwork samples.²

References

1. Lim, H.S., Park, J.J., Ko, K., *et al.* Syntheses of S1P analogues and their interaction with EDG/S1P receptors. *Bioor. Med. Chem. Lett.* **14(10)**, 2499-2503 (2004).
2. Aljohani, A.J., Edwards, G., Guerra, Y., *et al.* Human trabecular meshwork sphingolipid and ceramide profiles and potential latent fungal commensalism. *Invest. Ophthalmol. Vis. Sci.* **55(6)**, 3413-3422 (2014).

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, 05/01/2024

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