

PRODUCT INFORMATION



Sphingosine-1-phosphate (d18:1(14Z))

Item No. 27020

Formal Name: (2S,3R,Z)-2-amino-3-hydroxyoctadec-14-en-1-yl dihydrogen phosphate

Synonym: S1P (d18:1(14Z))

MF: C₁₈H₃₈NO₅P

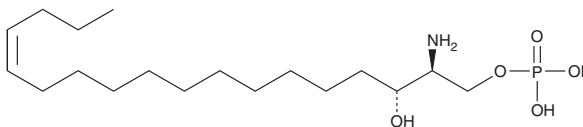
FW: 379.5

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

Sphingosine-1-phosphate (d18:1(14Z)) is supplied as a crystalline solid. Aqueous solutions of sphingosine-1-phosphate (d18:1(14Z)) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of sphingosine-1-phosphate (d18:1(14Z)) in 0.3 M sodium hydroxide is approximately 4 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Sphingosine-1-phosphate (d18:1(14Z)) is an atypical sphingolipid that contains a *cis* double bond at the 14-15 position rather than a *trans* double bond at the 4-5 position as in sphingosine-1-phosphate (d18:1) (Item No. 62570).

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, 02/16/2024

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