

PRODUCT INFORMATION



Sphingosine-1-phosphate (d18:1) Alkyne

Item No. 28965

Formal Name: (2S,3R,E)-2-amino-3-hydroxyoctadec-4-en-17-yn-1-yl dihydrogen phosphate

Synonyms: Click Tag™ Sphingosine-1-phosphate (d18:1) Alkyne, S1P (d18:1) Alkyne, Sphingosine-1-Phosphoric Acid Alkyne

MF: C₁₈H₃₄NO₅P

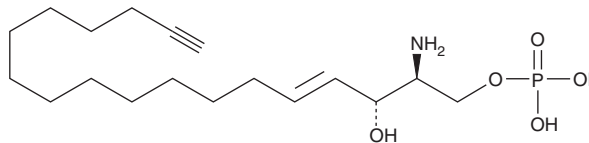
FW: 375.4

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

Sphingosine-1-phosphate (d18:1) alkyne is supplied as a crystalline solid. A stock solution may be made by dissolving the sphingosine-1-phosphate (d18:1) alkyne in the solvent of choice, which should be purged with an inert gas. Sphingosine-1-phosphate (d18:1) alkyne is soluble in 0.3 M NaOH at a concentration of approximately 4 mg/ml.

Description

Sphingosine-1-phosphate alkyne is an ω-alkyne form of sphingosine-1-phosphate (Item No. 62570). The terminal alkyne group can be used in click chemistry reactions to tag sphingosine-1-phosphate with fluorescent or biotinylated labels for analysis of its metabolism and biological activity.

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, 03/20/2024

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