

PRODUCT INFORMATION



Biotin Sphingosine

Item No. 39291

CAS Registry No.: 752987-57-8
Formal Name: (3a*S*,4*S*,6a*R*)-*N*-[(14*E*,16*R*,17*S*)-17-amino-16,18-dihydroxy-14-octadecen-1-yl]hexahydro-2-oxo-1*H*-thieno[3,4-*d*]imidazole-4-pentanamide

Synonym: Biotinyl Sphingosine

MF: C₂₈H₅₂N₄O₄S

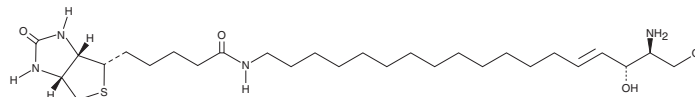
FW: 540.8

Purity: ≥95%

Supplied as: A 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

Biotin sphingosine is supplied as a solid. A stock solution may be made by dissolving the biotin sphingosine in the solvent of choice, which should be purged with an inert gas. Biotin sphingosine is soluble in methanol.

Description

Biotin sphingosine is a biotinylated derivative of sphingosine (d18:1) (Item No. 10007907).¹ It has been used in assays to detect the phosphorylation activity of sphingosine kinase 1 (SPHK1) and SPHK2 in cell lysates.

Reference

1. Roberts, J.L., Moretti, P.A.B., Darrow, A.L., *et al.* An assay for sphingosine kinase activity using biotinylated sphingosine and streptavidin-coated membranes. *Anal. Biochem.* **331(1)**, 122-129 (2004).

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, 07/25/2023

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