

PRODUCT INFORMATION



C6 Biotin dihydro Ceramide (d18:0/6:0)

Item No. 28949

Formal Name: N-((2S,3R)-1,3-dihydroxyoctadecan-2-yl)-6-(5-((3aS,4S,6aR)-2-oxohexahydro-1H-thieno[3,4-d]imidazol-4-yl)pentanamido)hexanamide

Synonyms: Biotin Ceramide (d18:0/6:0), Biotin Cer(d18:0/6:0), N-hexanoyl-biotin-D-erythro-Dihydrosphingosine

MF: C₃₄H₆₄N₄O₅S

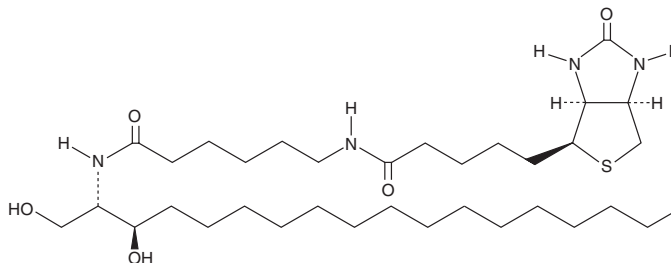
FW: 641.0

Purity: ≥98%

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

C6 Biotin dihydro ceramide (d18:0/6:0) is supplied as a solid. A stock solution may be made by dissolving the C6 biotin dihydro ceramide (d18:0/6:0) in the solvent of choice, which should be purged with an inert gas. C6 Biotin dihydro ceramide (d18:0/6:0) is soluble in a 2:1 solution of chloroform:methanol.

Description

C6 Biotin dihydro ceramide is an affinity probe that allows C6 dihydro ceramide (Item No. 24368) to be detected or immobilized through interaction with the biotin ligand.

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, 12/06/2022

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