

PRODUCT INFORMATION



C6 Phytoceramide (t18:0/6:0)

Item No. 24438

CAS Registry No.: 249728-94-7

Formal Name: N-[(1S,2S,3R)-2,3-dihydroxy-1-(hydroxymethyl)heptadecyl]-hexanamide

Synonyms: Cer(t18:0/6:0), Ceramide (t18:0/6:0), C6 Phytosphingosine (t18:0/6:0), N-hexanoyl Phytosphingosine,

MF: $C_{24}H_{49}NO_4$

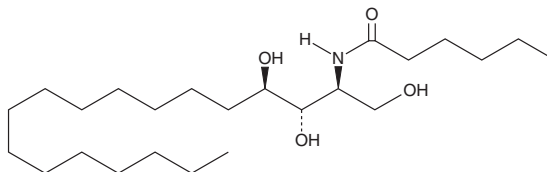
FW: 415.7

Purity: $\geq 98\%$

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 Phytoceramide (t18:0/6:0) is supplied as a solid. A stock solution may be made by dissolving the C6 phytoceramide (t18:0/6:0) in the solvent of choice. C6 Phytoceramide (t18:0/6:0) is soluble in ethanol and methanol and a 1:1 (warm) solution of chloroform:methanol.

Description

C6 Phytoceramide is a cell-permeable analog of phytoceramide (Item No. 20217). It is composed of a phytoceramide backbone amine-linked to a C6 fatty acid chain. C6 Phytoceramide induces the expression of differentiation-associated genes and promotes differentiation in keratinocytes *in vitro*.¹

Reference

1. Paragh, G., Schling, P., Ugocsai, P., et al. Novel sphingolipid derivatives promote keratinocyte differentiation. *Exp. Dermatol.* **17(12)**, 1004-1016 (2008).

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