

PRODUCT INFORMATION



C23 Phytoceramide (t18:0/23:0)

Item No. 9004471

CAS Registry No.: 844866-89-3

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

Synonyms: C23 Phytosphingosine (t18:0/23:0), Ceramide (t18:0/23:0), Cer(t18:0/23:0), N-Tricosanoyl Phytosphingosine

MF: $C_{41}H_{83}NO_4$

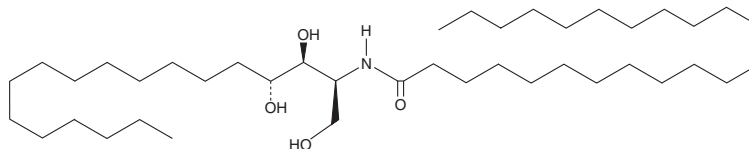
FW: 654.1

Purity: $\geq 98\%$

Supplied as: A solid

Storage: -20°C

Stability: ≥ 2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

C23 Phytoceramide (t18:0/23:0) is a long-chain sphingolipid that has been found in *C. sinensis*.¹

Reference

1. Valsecchi, M., Mauri, L., Casellato, R., *et al.* Ceramides as possible nutraceutical compounds: Characterization of the ceramides of the Moro blood orange (*Citrus sinensis*). *J. Agric. Food Chem.* **60**(40), 10103-10110 (2012).

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, 08/12/2024

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