

PRODUCT INFORMATION



C22 Phytoceramide-d₃ (t18:0/22:0-d₃)

Item No. 9003465

CAS Registry No.: 2011762-87-9

Formal Name: N-[(1S)-2S,3R-dihydroxy-1-(hydroxymethyl)heptadecyl]-docosanamide-d₃

Synonyms: Ceramide (t18:0/22:0-d₃),
Cer(t18:0/22:0-d₃),
C22 Phytosphingosine-d₃ (t18:0/22:0-d₃),
N-Docosanoyl Phytosphingosine-d₃

MF: C₄₀H₇₈D₃NO₄

FW: 643.1

Chemical Purity: ≥98% (C22 Phytoceramide (t18:0/22:0))

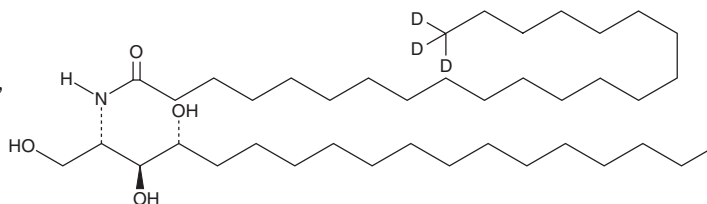
Deuterium

Incorporation: ≥99% deuterated forms (d₁-d₃); ≤1% d₀

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

C22 Phytoceramide-d₃ is intended for use as an internal standard for the quantification of C22 phytoceramide (Item No. 9003464) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated *versus* unlabeled).

Description

C22 Phytoceramide is a sphingolipid that has been found in Moro blood orange (*C. sinensis*) peels.¹ It has also been found in primary human astrocytes.²

References

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).
2. Dasgupta, S., Kong, J., and Bieberich, E. Phytoceramide in vertebrate tissues: One step chromatography separation for molecular characterization of ceramide species. *PLoS One* **8**(11), e80841 (2013).

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, 12/13/2022

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