

# PRODUCT INFORMATION



## C22 Phytoceramide (t18:0/22:0)

Item No. 9003464

**CAS Registry No.:** 164576-03-8  
**Formal Name:** N-[(1S,2S,3R)-2,3-dihydroxy-1-(hydroxymethyl)heptadecyl]-docosanamide

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

**MF:** C<sub>40</sub>H<sub>81</sub>NO<sub>4</sub>

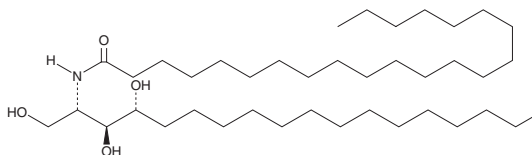
**FW:** 640.1

**Purity:** ≥98%

**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.

### Description

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

### 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, 11/30/2022

#### 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