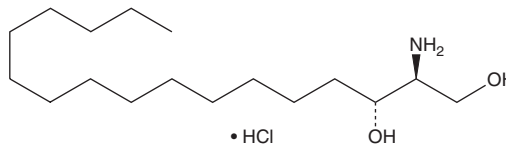


PRODUCT INFORMATION

Sphinganine (d17:0) (hydrochloride)

Item No. 37989

Formal Name: 2S-amino-1,3R-heptadecanediol, monohydrochloride
Synonyms: D-erythro-C17-Dihydrosphingosine, D-erythro-Sphinganine C-17
MF: $C_{17}H_{37}NO_2 \cdot HCl$
FW: 323.9
Purity: $\geq 95\%$
Supplied as: A crystalline solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



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

Laboratory Procedures

Sphinganine (d17:0) (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the sphinganine (d17:0) (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Sphinganine (d17:0) (hydrochloride) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of sphinganine (d17:0) (hydrochloride) in these solvents is approximately 2 and 10 mg/ml, respectively.

Description

Sphinganine (d17:0) is a synthetic bioactive sphingolipid.¹ It is active against *C. glabrata* and *C. albicans* with a minimum fungicidal concentration (MFC) value of 0.5 $\mu g/ml$ for both. Sphinganine (d17:0) has commonly been used as an internal standard for the detection of sphingolipids in chromatographic and mass spectrometry applications.^{2,3}

References

1. Thevissen, K., Hillaert, U., Meert, E.M., *et al.* Fungicidal activity of truncated analogues of dihydrosphingosine. *Bioorg. Med. Chem. Lett.* **18(13)**, 3728-3730 (2008).
2. Spassieva, S., Bielawski, J., Anelli, V., *et al.* Combination of C_{17} sphingoid base homologues and mass spectrometry analysis as a new approach to study sphingolipid metabolism. *Methods Enzymol.* **434**, 233-241 (2007).
3. Qu, F., Wu, C.S., Hou, J.F., *et al.* Sphingolipids as new biomarkers for assessment of delayed-type hypersensitivity and response to triptolide. *PLoS One* **7(12)**, e52454 (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, 03/27/2024

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