

PRODUCT INFORMATION



DL-erythro Sphinganine (d20:0)

Item No. 24378

Formal Name: 2R-aminoicosane-1,3S-diol,
2R-monoaminoicosane-1,3S-diol

Synonyms: DL-erythro Dihydrosphingosine (d20:0),
C20 DL-erythro Sphinganine

MF: C₄₀H₈₆N₂O₄

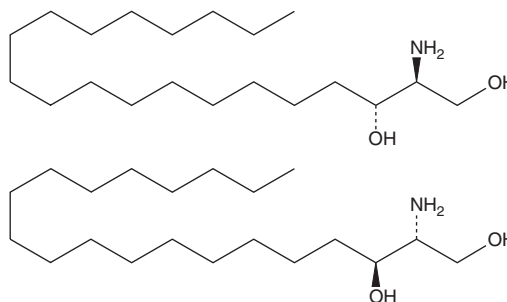
FW: 659.1

Purity: ≥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

DL-erythro Sphinganine (d20:0) is supplied as a solid. A stock solution may be made by dissolving the DL-erythro sphinganine (d20:0) in the solvent of choice. DL-erythro Sphinganine (d20:0) is soluble in a 5:1 solution of chloroform:methanol. DL-erythro Sphinganine (d20:0) is also soluble in ethanol (warmed).

Description

DL-erythro Sphinganine (d20:0) is a mixture of naturally occurring sphinganine (d20:0) (Item No. 22509) and synthetic L-erythro sphinganine (d20:0).

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, 10/31/2022

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