

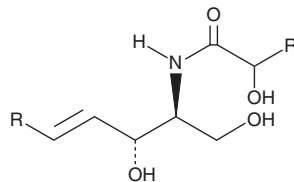
PRODUCT INFORMATION



Ceramides Mixture

Item No. 22853

CAS Registry No.: 100403-19-8
Purity: ≥98% (mixture of ceramides)
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

Ceramides mixture is supplied as a crystalline solid. A stock solution may be made by dissolving the ceramides mixture in the solvent of choice, which should be purged with an inert gas. Ceramides mixture is soluble in the organic solvent chloroform at a concentration of approximately 10 mg/ml.

Description

The ceramides mixture consists of hydroxy and non-hydroxy fatty acid-containing ceramides. Ceramides are generated from sphingomyelin through activation of sphingomyelinases or through the *de novo* synthesis pathway, which requires the coordinated action of serine palmitoyl transferase and ceramide synthase. They have been shown to mediate antiproliferative responses such as apoptosis, growth arrest, differentiation, and senescence.¹

Reference

1. Ogretmen, B., Pettus, B.J., Rossi, M.J., *et al.* Biochemical mechanisms of the generation of endogenous long chain ceramide in response to exogenous short chain ceramide in the A547 human lung adenocarcinoma cell line. *J. Biol. Chem.* **277(15)**, 12960-12969 (2002).

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, 10/27/2022

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