

PRODUCT INFORMATION



Sphingomyelinase

Item No. 10010076

Overview

Contents: This vial contains a lyophilized powder of sphingomyelinase from *Bacillus cereus*.
Storage: 4°C
Stability: ≥2 years

Laboratory Procedures

Prior to use in the Sphingomyelin Colorimetric Assay Kit (Item No. 10009928), reconstitute the Sphingomyelinase with 1 ml of diluted SM Buffer (10X) (Item No. 10010072). Store the reconstituted enzyme mixture on ice until ready to use. Unused enzyme mixture should be frozen at -20°C and will be stable for one month.

Precaution

This product may not perform as described if used with other assay reagents or protocols.

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, 01/19/2026

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