

PRODUCT INFORMATION



SM Enzyme Mixture

Item No. 10010074

Overview

Contents: This vial contains a lyophilized powder consisting of a mixture of choline oxidase and horseradish peroxidase.
Storage: 4°C
Stability: ≥1 year

Laboratory Procedures

Prior to used in the Sphingomyelin Colorimetric Assay Kit (Item No. 10009928), reconstitute the Enzyme Mixture contents with 1 ml of diluted SM Buffer (10X) (Item No. 10010072). Store the reconstituted Enzyme Mixture on ice until ready to use.

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, 09/28/2023

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