

PRODUCT INFORMATION

PC-Specific Phospholipase D

Item No. 10010069

Overview

Contents: This vial contains 100 ml of a phosphatidylcholine-specific phospholipase D (PLD) solution.
Storage: -20°C
Stability: ≥1 year

Laboratory Procedures

The PC-specific PLD is ready to use as supplied in the Phosphatidylcholine Colorimetric Assay Kit (Item No. 10009926).

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, 09/24/2023

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