

PRODUCT INFORMATION



PC Buffer (10X)

Item No. 10010066

Overview

Contents: This vial contains 5 ml of 500 mM Tris-HCl, pH 8.0, containing 6.6 mM calcium chloride.
Storage: -20°C
Stability: ≥1 year

Laboratory Procedures

Prior to use in the Phosphatidylcholine Colorimetric Assay Kit (Item No. 10009926), dilute 3 ml of Buffer concentrate with 27 ml of HPLC-grade water. This final Buffer (50 mM Tris-HCl, pH 8.0, containing 0.66 mM CaCl₂) should be used for reconstituting the PC Color Detector (Item No. 10010067), PC Enzyme Mixture (Item No. 10010068), and diluting the Phosphatidylcholine Standard (Item No. 10010070). When stored at 4°C, the diluted Buffer is stable for at least six months.

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, 09/29/2023

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