

PRODUCT INFORMATION



Glucose Colorimetric Enzyme Mixture

Item No. 10010100

Overview

Contents: This vial contains a lyophilized enzyme mixture.
Storage: -20°C
Stability: ≥1 year

Laboratory Procedures

Prior to performing the Glucose Colorimetric Assay Kit (Item No. 10009582), reconstitute the lyophilized enzyme mixture with 6 ml of diluted Sodium Phosphate Assay Buffer (Item No. 700003) and mix well. The reconstituted solution is stable for one hour when stored on ice. One vial of enzyme mixture is sufficient to evaluate 60 wells.

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, 09/22/2023

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