

# Product Information



## Glycolysis Assay Enzyme Mixture

Item No. 600452

**Contents:** This vial contains a lyophilized powder of enzymes.

**Stability:** ≥1 year at -20°C

### Laboratory Procedures

Prior to use in the Glycolysis Assay Kit (Item No. 600450), reconstitute the contents of the vial with 150 µl of Cell-Based Assay Buffer (Item No. 10009322) and put the vial on ice. The reconstituted mixture is stable for two hours at 4°C. If the reconstituted enzyme will not be used all at once, it is recommended that aliquots be frozen at -80°C. Freezing and thawing of the solution should be limited to a single time.

To prepare 12 ml of reaction solution enough for a 96-well plate assay, add 120 µl of this reconstitute enzyme solution along with 120 µl of 100 X Glycolysis Assay Substrate (Item No. 600451) and 120 µl of 100X Glycolysis Assay Cofactor (Item No. 600454) to 11.64 ml of Cell-Based Assay Buffer (Item No. 10009322).

### CAUTIONS

For laboratory research use only. This reagent is not for human or veterinary use. This product has been tested and formulated to work with the Glycolysis Cell-Based Assay Kit and Oxygen Consumption/Glycolysis Dual Assay Kit from Cayman (Item Nos. 600450 and 601060, respectively). This product may not perform as described if used with other assay reagents or protocols.

### Related Products

For a list of related products please visit: [www.caymanchem.com/catalog/600452](http://www.caymanchem.com/catalog/600452)

### Cayman Chemical

#### Mailing address

1180 E. Ellsworth Road  
Ann Arbor, MI  
48108 USA

#### Phone

(800) 364-9897  
(734) 971-3335

#### Fax

(734) 971-3640

#### E-Mail

[custserv@caymanchem.com](mailto:custserv@caymanchem.com)

#### Web

[www.caymanchem.com](http://www.caymanchem.com)

**WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.**

#### SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. 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

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery**.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

**For further details, please refer to our Warranty and Limitation of Remedy located on our website and in our catalog.**

Copyright Cayman Chemical Company, 02/25/2014