

# Product Information



## Cell-Based Assay Glutathione Standard

Item No. 600362

**Contents:** This vial contains 3 mg of lyophilized glutathione (GSH).

**Stability:**  $\geq 1$  year at  $-20^{\circ}\text{C}$

### Laboratory Procedures

For standard preparation, dissolve the content of the Cell-Based Assay Glutathione Standard in 1 ml of Cell-Based Assay Buffer (prepared from Cell-Based Assay Buffer Tablet, Item No. 10009322). The concentration of this GSH solution will be 10 mM. This GSH Standard is stable for four weeks at  $4^{\circ}\text{C}$ .

For standard preparation as a positive control in a standard curve, dilute the contents of the vial 1:100 in Assay Buffer. Load 90  $\mu\text{l}$  of this control into corresponding wells in the assay plate.

To run a GSH standard curve, obtain eight clean test tubes and label them #1 to # 8. Add 975  $\mu\text{l}$  of Assay Buffer into tube #1 and 250  $\mu\text{l}$  into tubes #2-#8. Transfer 25  $\mu\text{l}$  of the GSH Standard from the 10 mM GSH Standard Solution into tube #1 and mix thoroughly. The concentration of this standard, the first point on the standard curve, is 250  $\mu\text{M}$ . Serially dilute the standard by removing 250  $\mu\text{l}$  from tube #1 and placing it into tube #2; mix thoroughly. Next remove 250  $\mu\text{l}$  from tube #2 and place it into tube #3; mix thoroughly. Repeat it for tubes #4 and #7. Do not add any standard to tube #6. This tube will be your blank.

### CAUTIONS

For laboratory research use only. This reagent is not for human or veterinary use. This product has been tested and formulated to work exclusively with the Glutathione Cell-Based Detection Kit (Blue Fluorescence) from Cayman (Item No. 600360). This product may not perform as described if used with other assay reagents or protocols.

### Related Product

Glutathione Cell-Based Detection Kit (Blue Fluorescence) - Item No. 600360

## 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

### Web

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.**

#### MATERIAL 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 Material 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, 01/11/2011