

# PRODUCT INFORMATION

## GSH Co-Factor Mixture

Item No. 703016



### Overview

---

**Contents:** This vial contains a lyophilized powder of NADP<sup>+</sup> and glucose-6-phosphate.  
**Storage:** 4°C  
**Stability:** ≥1 years

### Laboratory Procedures

---

GSH Co-Factor Mixture has been tested and formulated to work exclusively with Cayman's Glutathione Assay Kit (Item No. 703002). Please visit [www.caymanchem.com/catalog/703002](http://www.caymanchem.com/catalog/703002) for the kit protocol, procedures, and product handling.

### 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, 12/02/2021

#### 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](http://WWW.CAYMANCHEM.COM)