

# PRODUCT INFORMATION

## Hemoglobin Detector (4X)

Item No. 700542



### Overview

---

**Contents:** This vial contains a solution of concentrated hemoglobin detector.  
**Storage:** 4°C  
**Stability:** ≥3 years

### Laboratory Procedures

---

Prior to use in the Hemoglobin Colorimetric Assay Kit (Item No. 700540), dilute the contents of the vial with 30 ml of HPLC-grade water. This is enough detector to assay one plate. When stored at 4°C, the diluted Hemoglobin Detector is stable for at least one month.

### 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, 11/21/2025

#### 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