

PRODUCT INFORMATION

Hemoglobin Sample Buffer (10X)

Item No. 700541



Overview

Contents: This vial contains a solution of 1 M potassium phosphate, pH 7.4, containing 10 mM EDTA.
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 90 ml of HPLC-grade water. This diluted Sample Buffer (100 mM potassium phosphate, pH 7.4, containing 1 mM EDTA) is used for dilution of samples. When stored at 4°C, the diluted buffer is stable for six months.

Precaution

For laboratory research use only. This reagent is not for human or veterinary use.

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, 11/20/2025

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