

PRODUCT INFORMATION

Potassium Phosphate Buffer (500 mM; pH 7.0)

Item No. 700621

Contents: The vial contains 5 ml of 500 mM potassium phosphate, pH 7.0.

Stability: ≥2 years at -20°C

Laboratory Procedures

Prior to use, dilute the contents of the vial with 45 ml of HPLC-grade water. This final Assay Buffer (50 mM potassium phosphate, pH 7.0) is stable for three months at 4°C.

Cautions

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, 12/10/2018

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