

PRODUCT INFORMATION



Uric Acid Standard

Item No. 700325

Overview

Contents: This vial contains a lyophilized powder of uric acid
Storage: -20°C
Stability: ≥1 year

Laboratory Procedures

Prior to use in the Uric Acid Assay Kit (Item No. 700320), add 1 ml of 1X Uric Acid Assay Buffer (Item No. 700321) to the vial and vortex. The concentration of the uric acid stock solution is 0.5 mM. It is ready to use to prepare the standard curve. The reconstituted standard is stable for four hours at room temperature. Please visit [www.caymanchem.com/catalog/kit number 700320](http://www.caymanchem.com/catalog/kit%20number%20700320) 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, 06/13/2017

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