

PRODUCT INFORMATION

Creatinine (serum) Color Reagent

Item No. 700461

Overview and Properties

Contents: The vial contains 12 ml of 1.2% picric acid.
Storage: Room temperature
Stability: ≥2 years

Procedures

The Color Reagent is ready to use as supplied in the Creatinine (serum) Colorimetric Assay Kit (Item No. 700460).

Cautions

For laboratory research use only. The Creatinine (serum) Color Reagent contains picric acid and thus it is recommended to take appropriate precautions when using the reagent (i.e., gloves, safety glasses, lab coat, etc.). It is harmful if swallowed and irritating to eyes, respiratory system, and skin. In case of contact with skin or eyes, rinse immediately with plenty of water for 15 minutes. This reagent is not for human or veterinary use. This product has been tested and formulated to work exclusively with the Creatinine (serum) Colorimetric Assay Kit from Cayman (Item No. 700460). 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, 10/06/2021

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