

PRODUCT INFORMATION

Protein Carbonyl Guanidine Hydrochloride

Item No. 10005848



Overview

Contents: This vial contains a solution of guanidine hydrochloride.
Storage: 4°C
Stability: ≥2 years

Laboratory Procedures

Protein Carbonyl Guanidine Hydrochloride has been tested and formulated to work exclusively with Cayman's Protein Carbonyl Colorimetric Assay Kit (Item No. 10005020). Please visit www.caymanchem.com/catalog/10005020 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, 04/12/2022

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