

# PRODUCT INFORMATION

## Xanthine Oxidase Sample Buffer (10X)

Item No. 10010973



### Overview

**Contents:** The vial contains 5 ml of 1 M Tris-HCl, pH 7.5.  
**Storage:** -20°C  
**Stability:** ≥2 years

### Laboratory Procedures

Prior to use in the Xanthine Oxidase Fluorometric Assay Kit (Item No. 10010895), dilute 3 ml of Sample Buffer concentrate with 27 ml of HPLC-grade water. This final Sample Buffer (100 mM Tris-HCl, pH 7.5) is used for the preparation of the Xanthine Oxidase Standard (Item No. 10010974). The diluted Sample Buffer is stable for six months at 4°C.

### Precaution

For laboratory research use only. This reagent is not for human or veterinary use. This product has been tested and formulated to work exclusively with the Xanthine Oxidase Fluorometric Assay Kit from Cayman (Item No. 10010895). 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, 06/18/2026

#### 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](http://WWW.CAYMANCHEM.COM)