

CERTIFICATE of ANALYSIS



Xanthine Oxidase (bovine)

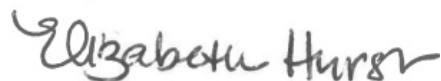
Item No. 38017 • Batch No. 0809371 • Expires 04MAY2027

Overview

Formulation:	2.3 M of ammonium sulfate with 10 mM sodium phosphate, pH 7.8, 1 mM EDTA and 1 mM sodium salicylate
Concentration:	(no concentration)
Storage/stability:	4°C/1yr
Unit definition:	One unit is defined as the amount of enzyme required to form 1 μ mol of uric acid per minute at room temperature in 50 mM Tris buffer, pH 8.0, using 0.1 mM hypoxanthine as a substrate
Activity:	1.35 U/ml

Reviewed and approved by: Beth Hurst

Approved date: 05MAY2026



WARNING

THIS PRODUCT IS FOR RESEARCH USE - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE. IT IS THE RESPONSIBILITY OF THE PURCHASER TO DETERMINE SUITABILITY FOR OTHER APPLICATIONS.

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, 10/12/2018

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