

PRODUCT INFORMATION



Xanthine Oxidase (bovine)

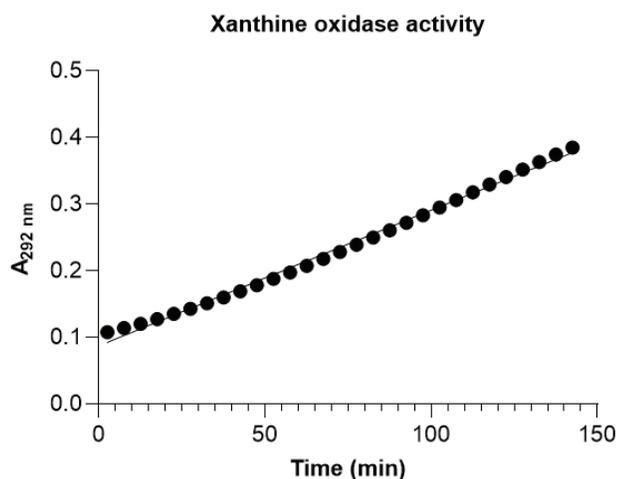
Item No. 38017

Overview and Properties

Synonyms:	Xanthine Dehydrogenase/Oxidase, Xanthine Oxidoreductase, XO, XOR
Source:	Raw bovine milk
Amino Acids:	1-1,332 (full length)
Molecular Weight:	146.79 kDa
Storage:	4°C (as supplied)
Stability:	≥6 months
Supplied in:	2.3 M Ammonium sulfate with 10 mM sodium phosphate, pH 7.8, 1 mM EDTA and 1 mM sodium salicylate
Activity:	<i>batch specific</i> U/ml
Unit Definition:	One unit is defined as the amount of enzyme required to formation of 1 μmol of uric acid per minute at room temperature in 50 mM Tris, pH 8.0, buffer using 0.1 mM hypoxanthine as a substrate.

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Image



Xanthine oxidase activity measured at 292 nm in the presence of 0.1 mM hypoxanthine and 0.1 mM DTPA at room temperature in 50 mM Tris, pH 8.0.

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, 07/31/2025

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

PRODUCT INFORMATION



Description

Xanthine oxidase (XO) is a complex metalloflavoprotein involved in purine catabolism, oxidative stress, and xenobiotic metabolism, among other processes.^{1,2} It exists as a homodimer and each monomer is composed of an N-terminal domain that contains two unequal iron-sulfur clusters, an intermediate domain that contains the FAD cofactor, and a C-terminal domain that contains the molybdopterin cofactor. XO is primarily expressed in the liver and intestines, has been found in blood, serum, and milk, and localizes to the cytosol, cell membrane, and peroxisomes.¹ It is produced by irreversible or reversible post-translational modification of xanthine dehydrogenase (XD) *via* limited proteolysis or oxidation of XD thiol groups, respectively.^{1,2} XO catalyzes the conversion of hypoxanthine to xanthine to uric acid and uses molecular oxygen as the electron acceptor, compared with XD that uses NAD⁺, and this results in the generation of hydrogen peroxide and superoxide, which can react with nitric oxide (NO) to produce peroxynitrite.¹⁻³ In hypoxic conditions, XO converts nitrates and nitrite to NO.⁴ Increased blood levels of XO's enzymatic product uric acid, hyperuricemia, is associated with several diseases, including gout, cardiovascular disease, ischemia-reperfusion injury, obesity, and diabetes.^{1,2,5,6,7} XO has commonly been used in coupled enzyme assays to measure superoxide dismutase (SOD) activity.^{8,9} Cayman's Xanthine Oxidase (bovine) protein can be used for enzyme activity assays.

References

1. Battelli, M.G., Bolognesi, A., and Polito, L. Pathophysiology of circulating xanthine oxidoreductase: New emerging roles for a multi-tasking enzyme. *Biochim. Biophys. Acta.* **1842(9)**, 1502-1517 (2014).
2. Battelli, M.G., Bortolotti, M., Bolognesi, A., *et al.* Pro-aging effects of xanthine oxidoreductase products. *Antioxidants (Basel)* **9(9)**, 839 (2020).
3. Lee, C., Liu, X., and Zweier, J.L. Regulation of xanthine oxidase by nitric oxide and peroxynitrite. *J. Biol. Chem.* **275(13)**, 9369-9376 (2000).
4. Millar, T.M., Stevens, C.R., Benjamin, N., *et al.* Xanthine oxidoreductase catalyses the reduction of nitrates and nitrite to nitric oxide under hypoxic conditions. *FEBS Lett.* **427**, 225-228 (1998).
5. Garcia-Gil, M., Camici, M., Allegrini, S., *et al.* Emerging role of purine metabolizing enzymes in brain function and tumors. *Int. J. Mol. Sci.* **19(11)**, 3598 (2018).
6. Dawson, J., and Walters, M. Uric acid and xanthine oxidase: Future therapeutic targets in the prevention of cardiovascular disease? *Br. J. Clin. Pharmacol.* **62(6)**, 633-644 (2006).
7. Glantzounis, G.K., Tsimoyiannis, E.C., Kappas, A.M., *et al.* Uric acid and oxidative stress. *Curr. Pharm. Des.* **11(32)**, 4145-4151 (2005).
8. Sun, Y., Oberley, L.W., and Li, Y. A simple method for clinical assay of superoxide dismutase. *Clin. Chem.* **34(3)**, 497-500 (1988).
9. Okado-Matsumoto, A. and Fridovich, I. Subcellular distribution of superoxide dismutases (SOD) in rat liver. Cu, Zn-SOD in mitochondria. *J. Biol. Chem.* **276(46)**, 38388-38393 (2001).

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