

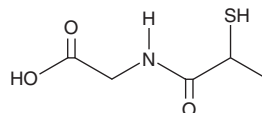
PRODUCT INFORMATION



Tiopronin

Item No. 23720

CAS Registry No.: 1953-02-2
Formal Name: N-(2-mercapto-1-oxopropyl)-glycine
Synonyms: BRN 1859822, (±)-Tiopronin
MF: C₅H₉NO₃S
FW: 163.2
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Tiopronin is supplied as a solid. A stock solution may be made by dissolving the tiopronin in the solvent of choice, which should be purged with an inert gas. Tiopronin is slightly soluble in DMSO and methanol.

Description

Tiopronin is an antioxidant that has diverse biological activities.¹⁻⁵ It reduces free radical production by murine macrophages and granulocytes *in vitro* in a dose-dependent manner.¹ Tiopronin induces expression of hypoxia-inducible factor 1α (HIF-1α) and increases VEGF secretion in human colon carcinoma cells.² Rectal administration (500 μL of a 10 mM solution) reduces myeloperoxidase activity and reduces pro-inflammatory cytokine production in the colon in a rat model of colitis. Tiopronin (80-320 mg/kg per day) reduces the incidence of cleft palate in fetuses born to female mice orally exposed to teratogenic methylmercury chloride.³ Tiopronin (100 mg/kg) reduces heme oxygenase 1 mRNA expression, lipid peroxidation, and transverse aortic constriction in a mouse model of cardiac hypertrophy.⁴ Tiopronin (20 mg/kg) is hepatoprotective, increasing activity of the antioxidant enzymes superoxide dismutase and glutathione peroxidase and reversing hepatocyte degeneration in a rat model of high-fat diet-induced non-alcoholic steatohepatitis.⁵ Formulations containing tiopronin have been used in the treatment of cystinuria, an inborn error of metabolism characterized by kidney stones and excessive urinary excretion of cysteine, arginine, ornithine, and lysine.

References

1. Rommain, M., Delecoeuillerie, G., Paul, J.L., *et al.* Inhibitory action of tiopronin on free radicals. *Rev. Rhum. Mal. Osteoartic.* **56**(5 Pt 2), 34-37 (1989).
2. Yum, S., Park, H., Hong, S., *et al.* N-(2-mercaptopropionyl)-glycine, a diffusible antioxidant, activates HIF-1 by inhibiting HIF prolyl hydroxylase-2: Implication in amelioration of rat colitis by the antioxidant. *Biochem. Biophys. Res. Commun.* **443**(3), 1008-1013 (2014).
3. Fujimoto, T., Fuyuta, M., Kiyofuji, E., *et al.* Prevention by tiopronin (2-mercaptopropionyl glycine) of methylmercuric chloride-induced teratogenic and fetotoxic effects in mice. *Teratology* **20**(2), 297-301 (1979).
4. Date, M.O., Morita, T., Yamashita, N., *et al.* The antioxidant N-2-mercaptopropionyl glycine attenuates left ventricular hypertrophy in in vivo murine pressure-overload model. *J. Am. Coll. Cardiol.* **39**(5), 907-912 (2002).
5. Wang, J.Q., Zou, Y.H., Huang, C., *et al.* Protective effects of tiopronin against high fat diet-induced non-alcoholic steatohepatitis in rats. *Acta. Pharmacol. Sin.* **33**(6), 791-797 (2012).

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, 06/30/2023

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