

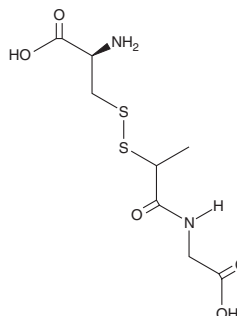
PRODUCT INFORMATION



Tiopronin-Cysteine Disulfide

Item No. 39020

CAS Registry No.: 77591-18-5
Formal Name: 3-[[2-[(carboxymethyl)amino]-1-methyl-2-oxoethyl]dithio]-L-alanine
Synonym: α-Mercaptopropionylglycine-Cysteine Disulfide
MF: C₈H₁₄N₂O₅S₂
FW: 282.3
Purity: ≥90%
Supplied as: A crystalline 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-cysteine disulfide is supplied as a crystalline solid. A stock solution may be made by dissolving the tiopronin-cysteine disulfide in water. We do not recommend storing the aqueous solution for more than one day.

Description

Tiopronin-cysteine disulfide is a metabolite of the antioxidant and anticystinuric agent tiopronin (Item No. 23720).¹

Reference

1. Backer, E.T., van den Ende, A., and Geus, W.P. Determination of urinary cystine for monitoring anticystinuric therapy. *Clin. Chim. Acta* **165(2-3)**, 351-358 (1987).

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, 06/22/2026

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