

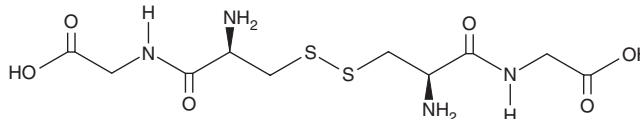
PRODUCT INFORMATION



Cystinyl-bis-glycine

Item No. 30494

CAS Registry No.: 7729-20-6
Formal Name: L-cysteinyl-glycine, bimol. (1→1')-disulfide
Synonyms: (Cys-Gly)₂, (H-Cys-Gly-OH)₂, NSC 333711
MF: C₁₀H₁₈N₄O₆S₂
FW: 354.4
Purity: ≥95%
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

Cystinyl-bis-glycine is supplied as a crystalline solid. Aqueous solutions of cystinyl-bis-glycine can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of cystinyl-bis-glycine in PBS, pH 7.2, is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Cystinyl-bis-glycine is a dipeptide intermediate in the synthesis of oxidized glutathione that is composed of two cysteinylglycine peptides linked by a disulfide bond.¹ It is formed *via* nonenzymatic oxidation of cysteinylglycine, a glutathione catabolite produced by γ -glutamyl transpeptidase.^{1,2}

References

1. Griffith, O.W. and Tate, S.S. The apparent glutathione oxidase activity of γ -glutamyl transpeptidase. Chemical mechanism. *J. Biol. Chem.* **255**(11), 5011-5014 (1980).
2. Kozak, E.M. and Tate, S.S. Glutathione-degrading enzymes of microvillus membranes. *J. Biol. Chem.* **257**(11), 6322-6327 (1982).

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, 12/12/2022

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