

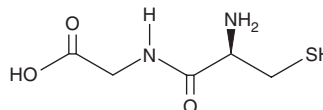
PRODUCT INFORMATION



Cysteinylglycine

Item No. 43262

CAS Registry No.: 19246-18-5
Formal Name: L-cysteinyl-glycine
Synonym: Cys-Gly
MF: $C_5H_{10}N_2O_3S$
FW: 178.2
Purity: $\geq 85\%$
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

Cysteinylglycine is supplied as a solid. A stock solution may be made by dissolving the cysteinylglycine in the solvent of choice, which should be purged with an inert gas. Cysteinylglycine is slightly soluble (0.1-1 mg/ml) in ethanol and DMSO.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of cysteinylglycine can be prepared by directly dissolving the solid in aqueous buffers. Cysteinylglycine is (soluble ≥ 10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

Cysteinylglycine is an endogenous dipeptide intermediate in glutathione (GSH) catabolism.¹ It is formed from extracellular GSH via hydrolysis or transpeptidation mediated by γ -glutamyl transpeptidase (GGT) and is a cellular source of cysteine and glycine for the formation of GSH.^{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. Zeevalk, G.D., Razmpour, R., and Bernard, L.P. Glutathione and Parkinson's disease: Is this the elephant in the room? *Biomed. Pharmacother.* **62**(4), 236-249 (2008).

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, 05/07/2025

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