

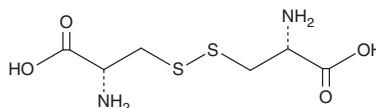
PRODUCT INFORMATION



L-Cystine

Item No. 31727

CAS Registry No.: 56-89-3
Synonyms: (-)-Cystine, NSC 13203
MF: C₆H₁₂N₂O₄S₂
FW: 240.3
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.

Description

L-Cystine is a dimeric form of cysteine that is formed by the covalent oxidative linkage of two cysteine residues.¹ It accumulates in the lysosomes of patients with cystinosis, an autosomal recessive lysosomal storage disorder, and is associated with renal Fanconi syndrome and loss of glomerular function.²

References

1. Appenzeller-Herzog, C. and Riemer, J. Techniques to monitor disulfide bond formation and the reduction potential of cysteine–cystine couples *in vitro* and *in vivo*. *Oxidative folding of proteins: Basic principles, cellular regulation and engineering*. Feige, M.J., editor, *The Royal Society of Chemistry* (2018).
2. Elmonem, M.A., Veys, K.R., Soliman, N.A., et al. Cystinosis: A review. *Orphanet J. Rare Dis.* **11**, 47 (2016).

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, 12/22/2022

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