

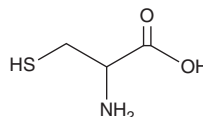
PRODUCT INFORMATION



DL-Cysteine

Item No. 41968

CAS Registry No.: 3374-22-9
Synonyms: (±)-Cysteine, (R,S)-Cysteine, NSC 63864
MF: C₃H₇NO₂S
FW: 121.2
Purity: ≥95%
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

DL-Cysteine is supplied as a solid. Aqueous solutions of DL-cysteine can be prepared by directly dissolving the solid in aqueous buffers. DL-Cysteine is soluble (≥10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

DL-Cysteine is a racemic mixture of D-cysteine and L-cysteine (Item No. 33309).¹ It prevents gastric ulceration induced by reserpine (Item No. 16474) and pyloric stenosis in rats when administered at a dose of 5% w/v.² Cryoprotectants containing DL-cysteine (5 mM) improve motility and DNA integrity in frozen and thawed epididymal spermatozoa from cats.³

References

1. Grishin, D.V., Zhdanov, D.D., Pokrovskaya, M.V., *et al.* D-amino acids in nature, agriculture and biomedicine. *All Life* **13(1)**, 11-22 (2020).
2. Salim, A.S. Protection against chronic gastric ulceration in the rat with sulfhydryl-containing agents. *J. Pharm. Sci.* **81(7)**, 698-700 (1992).
3. Thuwanut, P., Chatdarong, K., Techakumphu, M., *et al.* The effect of antioxidants on motility, viability, acrosome integrity and DNA integrity of frozen-thawed epididymal cat spermatozoa. *Theriogenology* **70(2)**, 233-240 (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, 12/17/2024

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