

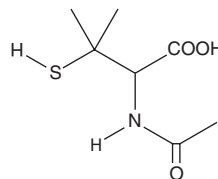
PRODUCT INFORMATION



N-Acetyl-DL-Penicillamine

Item No. 10010404

CAS Registry No.: 59-53-0
Formal Name: N-acetyl-3-mercapto-valine
Synonyms: NAPA, NSC 28039, NSC 92752
MF: C₇H₁₃NO₃S
FW: 191.2
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

N-Acetyl-DL-penicillamine is supplied as a crystalline solid. A stock solution may be made by dissolving the N-acetyl-DL-penicillamine in the solvent of choice, which should be purged with an inert gas. N-Acetyl-DL-penicillamine is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of N-acetyl-DL-penicillamine in these solvents is approximately 5 mg/ml in ethanol and 30 mg/ml in DMSO and DMF.

N-Acetyl-DL-penicillamine is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, N-acetyl-DL-penicillamine should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. N-Acetyl-DL-penicillamine has a solubility of approximately 0.12 mg/ml in a 1:7 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

N-Acetyl-DL-penicillamine is a chelating agent.¹⁻³ It inhibits the binding of methyl mercury to isolated human erythrocytes by 50% and removes 50% of methyl mercury ions from methyl mercury-loaded blood cells when used at a concentration of 1 mM.^{1,3} N-Acetyl-DL-penicillamine (3 mmol/kg per day, p.o.) reduces the biological half-life of mercury and decreases liver, kidney, brain, and blood mercury levels, as well as increases urinary excretion of mercury in a concentration-dependent manner, in mice when administered following injection of methyl mercuric chloride. It decreases mercuric chloride-induced mortality in mice when administered orally at a dose of 1.6 mmol/kg.² N-Acetyl-DL-penicillamine is also an analog of SNAP (Item No. 82250) that does not generate nitric oxide and has been used as a negative control in experiments using SNAP.^{4,5}

References

1. Aaseth, J. Mobilization of methyl mercury *in vivo* and *in vitro* using N-acetyl-DL-penicillamine and other complexing agents. *Acta Pharm. Toxicol. (Copenh.)* **39(3)**, 289-301 (1976).
2. Nielsen, J.B. and Andersen, O. Effect of four thiol-containing chelators on disposition of orally administered mercuric chloride. *Hum. Exp. Toxicol.* **10(6)**, 423-430 (1991).
3. Aaseth, J., Alexander, J., and Deverill, J. Evaluation of methyl mercury chelating agents using red blood cells and isolated hepatocytes. *Chem. Biol. Interact.* **36(3)**, 287-297 (1981).
4. Ogura, T., DeGeorge, G., Tatemichi, M., *et al.* Suppression of anti-microtubule agent-induced apoptosis by nitric oxide: Possible mechanism of a new drug resistance. *Jpn. J. Cancer Res.* **89(2)**, 199-205 (1998).
5. Takhampunya, R., Padmanabhan, R., and Ubol, S. Antiviral action of nitric oxide on dengue virus type 2 replication. *J. Gen. Virol.* **87(Pt. 10)**, 3003-3011 (2006).

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, 11/09/2022

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