

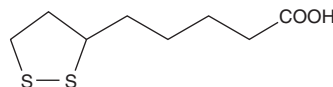
PRODUCT INFORMATION



DL- α -Lipoic Acid

Item No. 10005728

CAS Registry No.: 1077-28-7
Formal Name: 1,2-dithiolane-3-pentanoic acid
Synonym: Thiocctic Acid
MF: C₈H₁₄O₂S₂
FW: 206.3
Purity: \geq 95%
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

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

DL- α -lipoic acid is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, DL- α -lipoic acid should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. DL- α -lipoic acid has a solubility of approximately 0.25 mg/ml in a 1:8 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

DL- α -lipoic acid is a cyclic disulfide antioxidant that interconverts with its reduced dithiol form. It is an essential cofactor for decarboxylation reactions of the citric acid cycle, and acts as a general antioxidant.¹⁻³ DL- α -lipoic acid can act as a direct radical scavenger, as a cofactor to regenerate reduced glutathione, and as a metal chelator.

References

1. Biewenga, G.Ph., Haenen, G.R.M.M., and Bast, A. The pharmacology of the antioxidant lipoic acid. *General Pharmacology: The Vascular System* **29(3)**, 315-331 (1997).
2. Randle, P.J. Regulatory interactions between lipids and carbohydrates: The glucose fatty acid cycle after 35 years. *Diabetes Metab. Rev.* **14(4)**, 263-283 (1998).
3. Biewenga, G.Ph., Haenen, G.R.M.M., and Bast, A. The role of lipoic acid in the treatment of diabetic polyneuropathy*. *Drug Metab. Rev.* **29(4)**, 1025-1054 (1997).

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, 12/06/2022

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