

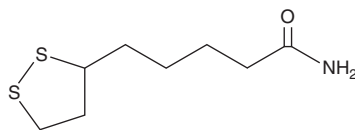
# PRODUCT INFORMATION



## Lipoamide

Item No. 35336

**CAS Registry No.:** 940-69-2  
**Formal Name:** 1,2-dithiolane-3-pentanamide  
**Synonyms:**  $\alpha$ -Lipoamide, DL- $\alpha$ -Lipoic Acid amide, NSC 90787, ( $\pm$ )-Thioctic Acid amide  
**MF:**  $C_8H_{15}NOS_2$   
**FW:** 205.3  
**Purity:**  $\geq 98\%$   
**Supplied as:** A solid  
**Storage:**  $-20^\circ\text{C}$   
**Stability:**  $\geq 4$  years



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

### Laboratory Procedures

Lipoamide is supplied as a solid. A stock solution may be made by dissolving the lipoamide in the solvent of choice, which should be purged with an inert gas. Lipoamide is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of lipoamide in these solvents is approximately 1, 2, and 5 mg/ml, respectively.

### Description

Lipoamide is an amide derivative of lipoic acid (Item No. 10005728).<sup>1</sup> It is an essential cofactor for enzymes in the citric acid cycle and enzymes involved in the catabolism of glycine and branched-chain amino acids (BCAAs).<sup>2</sup> Lipoamide is covalently attached to lysine residues in proteins via lipoylation, a post-translational modification that provides flexibility to enzyme complexes. It reduces hydrogen peroxide- or 6-OHDA-induced cytotoxicity in PC12 cells when used at concentrations of 50 and 100  $\mu\text{M}$  and increases the expression and protein levels of nuclear factor erythroid 2-related factor 2 (Nrf2) targets.<sup>1</sup> Lipoamide (100 mg/kg) also reverses mitochondrial dysfunction, increases the number of midbrain dopaminergic neurons, and improves fine motor activity in a rat model of Parkinson's disease induced by 6-OHDA.<sup>3</sup>

### References

1. Hou, Y., Li, X., Peng, S., *et al.* Lipoamide ameliorates oxidative stress via induction of Nrf2/ARE signaling pathway in PC12 cells. *J. Agric. Food Chem.* **67**(29), 8227-8234 (2019).
2. Rowland, E.A., Snowden, C.K., and Cristea, I.M. Protein lipoylation: An evolutionarily conserved metabolic regulator of health and disease. *Curr. Opin. Chem. Biol.* **42**, 76-85 (2018)
3. Zhou, B., Wen, M., Lin, X., *et al.* Alpha lipoamide ameliorates motor deficits and mitochondrial dynamics in the Parkinson's disease model induced by 6-hydroxydopamine. *Neurotox. Res.* **33**(4), 759-767 (2018).

#### 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, 11/23/2022

#### 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](http://WWW.CAYMANCHEM.COM)