

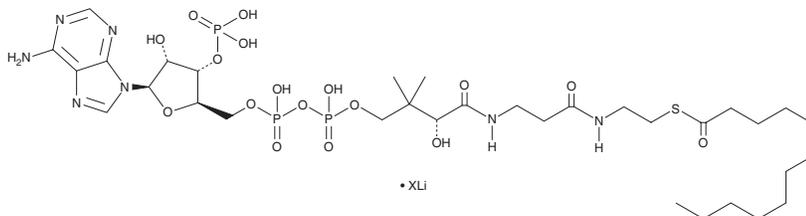
PRODUCT INFORMATION



Lauroyl-Coenzyme A (lithium salt)

Item No. 44593

CAS Registry No.: 190063-12-8
Formal Name: S-dodecanoate, coenzyme A, lithium salt
Synonyms: C12-CoA, Dodecanoyl-CoA, Dodecanoyl-Coenzyme A, Lauroyl-CoA
MF: C₃₃H₅₈N₇O₁₇P₃S • XLi
FW: 949.8
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Special Conditions: Light sensitive



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

Laboratory Procedures

Lauroyl-Coenzyme A (lauroyl-CoA) (lithium salt) is supplied as a solid. A stock solution may be made by dissolving the lauroyl-CoA (lithium salt) in the solvent of choice, which should be purged with an inert gas. Lauroyl-CoA (lithium salt) is sparingly soluble (1-10 mg/ml) in acetonitrile:water (1:1).

Description

Lauroyl-CoA is a medium-chain saturated fatty acid CoA, substrate for acyl-CoA thioesterases, and precursor to fatty acids.¹

Reference

1. Hunt, M.C., Ruitter, J., Mooyer, P., *et al.* Identification of fatty acid oxidation disorder patients with lowered acyl-CoA thioesterase activity in human skin fibroblasts. *Eur. J. Clin. Invest.* **35(1)**, 38-46 (2005).

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, 01/21/2026

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