

PRODUCT INFORMATION



Heptadecanoyl-Coenzyme A (lithium salt hydrate)

Item No. 25999

Formal Name: S-heptadecanoate coenzyme A, lithium salt hydrate

Synonym: Heptadecanoyl-CoA

MF: $C_{38}H_{68}N_7O_{17}P_3S \cdot XLi \cdot XH_2O$

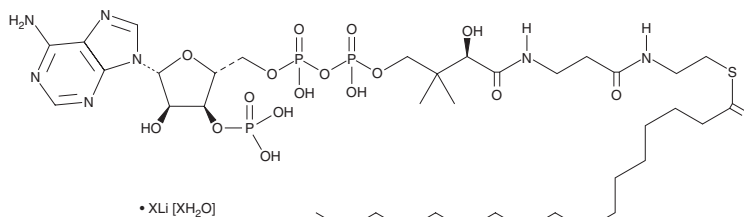
FW: 1,020.0

Purity: $\geq 95\%$

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

Heptadecanoyl-coenzyme A (heptadecanoyl-CoA) (lithium salt hydrate) is supplied as a solid. A stock solution may be made by dissolving the heptadecanoyl-CoA (lithium salt hydrate) in water. The solubility of heptadecanoyl-CoA (lithium salt hydrate) in water is approximately 20 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Heptadecanoyl-CoA is a long-chain saturated fatty acyl CoA. It has been used as an internal standard for the quantification of hepatic and mitochondrial fatty acyl CoAs.¹

Reference

1. Zhou, M., Xu, A., Lam, K.S.L., et al. Rosiglitazone promotes fatty acyl CoA accumulation and excessive glycogen storage in livers of mice without adiponectin. *J. Hepatol.* **53**(6), 1108-1116 (2010).

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, 12/20/2022

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