

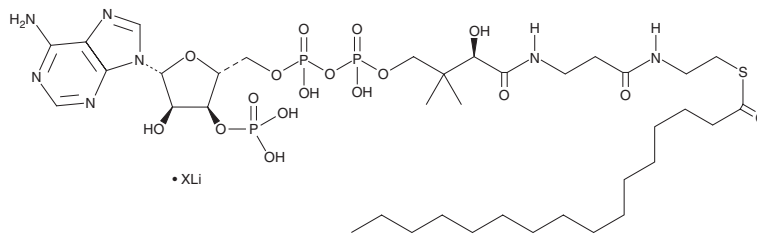
PRODUCT INFORMATION



Palmitoyl-Coenzyme A (lithium salt)

Item No. 45165

CAS Registry No.: 188174-64-3
Formal Name: S-hexadecanoate coenzyme A, lithium salt
Synonyms: Hexadecanoyl-CoA, Palmitoyl-CoA
MF: $C_{37}H_{66}N_7O_{17}P_3S \cdot XLi$
FW: 1,005.9
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

Palmitoyl-coenzyme A (lithium salt) is supplied as a solid. A stock solution may be made by dissolving the palmitoyl-coenzyme A (lithium salt) in water. Palmitoyl-coenzyme A (lithium salt) is soluble (≥ 10 mg/ml) in water. We do not recommend storing the aqueous solution for more than one day.

Description

Palmitoyl-coenzyme A (palmitoyl-CoA) is a derivative of CoA (Item Nos. 16147 | 21499) that contains the saturated fatty acid palmitic acid (Item Nos. 10006627 | 43561). It is a substrate for palmitoyltransferases during palmitoylation, a process that adds a palmitoyl group to proteins to anchor them to the membrane or target them to lipid rafts.¹ It is also a substrate in the *de novo* synthesis of sphingolipids.^{1,2} Palmitoyl-CoA binds to acyl-CoA-binding protein (ACBP; $K_d = 0.45$ pM) and undergoes β -oxidation in mitochondria.²

References

1. Guan, X. and Fierke, C. A. Understanding protein palmitoylation: Biological significance and enzymology. *Sci. China Chem.* **54**(12), 1888-1897 (2011).
2. Rasmussen, J. T., Faergeman, N. J., Kristiansen, K., *et al.* Acyl-CoA-binding protein (ACBP) can mediate intermembrane acyl-CoA transport and donate acyl-CoA for β -oxidation and glycerolipid synthesis. *Biochem. J.* **299**(Pt 1), 165-170 (1994)

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, 03/10/2026

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