

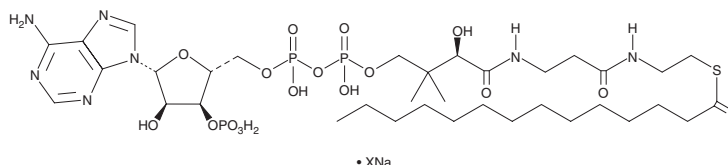
# PRODUCT INFORMATION



## Myristoyl-Coenzyme A (sodium salt)

Item No. 39777

**Formal Name:** S-tetradecanoate coenzyme A, sodium salt  
**Synonyms:** Myristoyl-CoA, Tetradecanoyl-CoA  
**MF:**  $C_{35}H_{62}N_7O_{17}P_3S \cdot XNa$   
**FW:** 977.9  
**Purity:**  $\geq 90\%$   
**Supplied as:** A solid  
**Storage:**  $-20^{\circ}C$   
**Stability:**  $\geq 4$  years



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

### Laboratory Procedures

Myristoyl-coenzyme A (myristoyl-CoA) (sodium salt) is supplied as a solid. A stock solution may be made by dissolving the myristoyl-CoA (sodium salt) in water. The solubility of myristoyl-CoA (sodium salt) in water is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Myristoyl-coenzyme A (myristoyl-CoA) is a derivative of CoA (Item No. 16147) that contains the saturated fatty acid myristic acid (Item No. 13351). It is a substrate for N-myristoyltransferase during myristoylation, a process that adds a myristoyl group to proteins either during translation to modify protein activity or post-translationally in apoptotic cells.<sup>1,2</sup> It is also a substrate in the *de novo* synthesis of phosphatidylinositol.<sup>3</sup>

### References

1. Bhatnagar, R.S., Fütterer, K., Waksman, G., *et al.* The structure of myristoyl-CoA: Protein N-myristoyltransferase. *Biochim. Biophys. Acta* **1441(2-3)**, 162-172 (1999).
2. Martin, D.D.O., Beauchamp, E., and Berthiaume, L.G. Post-translational myristoylation: Fat matters in cellular life and death. *Biochimie* **93(1)**, 18-31 (2011).
3. Darnell, J.C., Osterman, D.G., and Saltiel, A.R. Fatty acid remodelling of phosphatidylinositol under conditions of *de novo* synthesis in rat liver microsomes. *Biochim. Biophys. Acta* **1084(3)**, 279-291 (1991).

#### 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, 09/18/2023

#### 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