

PRODUCT INFORMATION



Pentanoyl-Coenzyme A (sodium salt)

Item No. 40460

Formal Name: coenzyme A, S-pentanoate, sodium salt

Synonyms: C5-CoA, Pentanoyl-CoA, Valeryl-CoA, Valeryl-Coenzyme A

MF: $C_{26}H_{44}N_7O_{17}P_3S \cdot XNa$

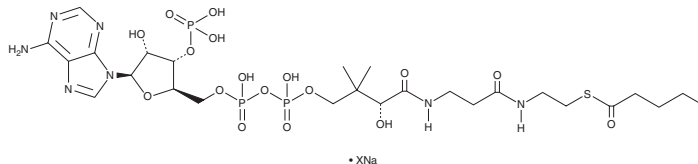
FW: 851.7

Purity: $\geq 95\%$

Supplied as: A solid

Storage: $-20^{\circ}C$

Stability: ≥ 4 years



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

Laboratory Procedures

Pentanoyl-coenzyme A (pentanoyl-CoA) (sodium salt) is supplied as a solid. A stock solution may be made by dissolving the pentanoyl-CoA (sodium salt) in water. We do not recommend storing the aqueous solution for more than one day.

Description

Pentanoyl-CoA is a derivative of CoA (Item Nos. 16147 | 21499 | 21722) that contains the short-chain saturated fatty acid pentanoic acid and a metabolite of 3-oxovalproyl-CoA.¹ It is formed from 3-oxovalproyl-CoA by mitochondrial thiolases. Hepatic levels of pentanoyl-CoA and propionyl-CoA are increased in an *Apc^{+/Min}-FCCC* mouse model of spontaneous colorectal cancer fed a low- or high-fat diet.²

References

1. Silva, M.F., Ruiter, J.P., Overmars, H., *et al.* Complete β -oxidation of valproate: Cleavage of 3-oxovalproyl-CoA by a mitochondrial 3-oxoacyl-CoA thiolase. *Biochem. J.* **362**(Pt 3), 755-760 (2002).
2. Dazard, J.-E., Sandlers, Y., Doerner, S.K., *et al.* Metabolomics of *Apc^{Min/+}* mice genetically susceptible to intestinal cancer. *BMC Syst. Biol.* **8:72**, (2014).

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, 03/13/2024

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