

PRODUCT INFORMATION



Octanoyl-Coenzyme A (sodium salt)

Item No. 27868

Formal Name: S-octanoate coenzyme A, sodium salt

Synonym: Octanoyl-CoA

MF: $C_{29}H_{50}N_7O_{17}P_3S \cdot XNa$

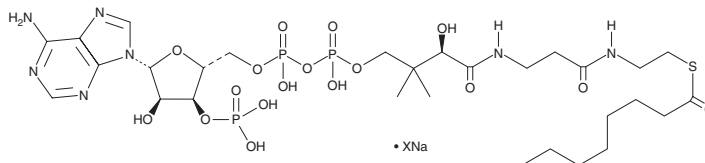
FW: 893.7

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

Octanoyl-coenzyme A (octanoyl-CoA) (sodium salt) is supplied as a solid. A stock solution may be made by dissolving the octanoyl-CoA (sodium salt) in the solvent of choice, which should be purged with an inert gas. Octanoyl-CoA (sodium salt) is soluble in organic solvents. Octanoyl-CoA (sodium salt) is slightly soluble in ethanol.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of octanoyl-CoA (sodium salt) can be prepared by directly dissolving the solid in aqueous buffers. The solubility of octanoyl-CoA (sodium salt) in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Octanoyl-CoA is a medium-chain acyl CoA and a metabolic intermediate in mitochondrial fatty acid β -oxidation.¹⁻³ Levels of octanoyl-CoA are increased in the liver of patients with Reye's syndrome and β -oxidation of octanoyl-CoA by medium-chain acyl CoA dehydrogenase (MCADH) is decreased in patients with MCADH deficiency (MCD).^{2,3} Octanoyl-CoA inhibits citrate synthase and glutamate dehydrogenase.¹

References

1. Lai, J.C., Liang, B.B., Jarvi, E.J., *et al.* Differential effects of fatty acyl coenzyme A derivatives on citrate synthase and glutamate dehydrogenase. *Res. Commun. Chem. Pathol. Pharmacol.* **82(3)**, 331-338 (1993).
2. Corkey, B.E., Hale, D.E., Glennon, M.V., *et al.* Relationship between unusual hepatic acyl coenzyme A profiles and the pathogenesis of Reye syndrome. *J. Clin. Invest.* **82(3)**, 782-788 (1988).
3. Amendt, B.A. and Rhead, W.J. Catalytic defect of medium-chain acyl-coenzyme A dehydrogenase deficiency. Lack of both cofactor responsiveness and biochemical heterogeneity in eight patients. *J. Clin. Invest.* **76(3)**, 963-969 (1985).

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, 10/13/2022

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