

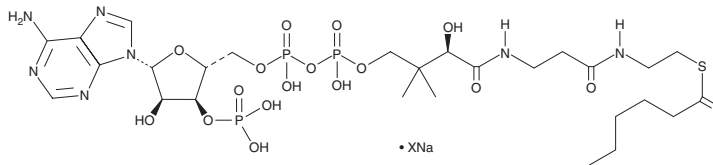
PRODUCT INFORMATION



Hexanoyl-Coenzyme A (sodium salt)

Item No. 27866

Formal Name: S-hexanoate coenzyme A, sodium salt
Synonyms: Caproyl-Coenzyme A, C6:0 CoA, Hexanoate-Coenzyme A
MF: $C_{27}H_{46}N_7O_{17}P_3S \cdot XNa$
FW: 865.7
Purity: $\geq 90\%$
UV/Vis.: λ_{max} : 258 nm
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years
Item Origin: Synthetic



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

Laboratory Procedures

Hexanoyl-coenzyme A (CoA) (sodium salt) is supplied as a solid. A stock solution may be made by dissolving the hexanoyl-CoA (sodium salt) in the solvent of choice, which should be purged with an inert gas. Hexanoyl-CoA (sodium salt) is slightly soluble in the organic solvent 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 hexanoyl-CoA (sodium salt) can be prepared by directly dissolving the solid in aqueous buffers. The solubility of hexanoyl-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

Hexanoyl-CoA is a medium-chain fatty acyl CoA. It is the preferred acyl donor substrate for ghrelin O-acyltransferase (GOAT) over octanoyl-CoA (Item No. 27868) in enzyme assays.¹ It is a precursor in the biosynthesis of olivetolic acid, a resorcinolic acid precursor in the synthesis of certain phytocannabinoids.²

References

- Ohgusu, H., Shirouzu, K., Nakamura, Y., *et al.* Ghrelin O-acyltransferase (GOAT) has a preference for n-hexanoyl-CoA over n-octanoyl-CoA as an acyl donor. *Biochem. Biophys. Res. Commun.* **386**(1), 153-158 (2009).
- Taura, F., Tanaka, S., Taguchi, C., *et al.* Characterization of olivetol synthase, a polyketide synthase putatively involved in cannabinoid biosynthetic pathway. *FEBS Lett.* **583**(12), 2061-2066 (2009).

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, 09/13/2023

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