

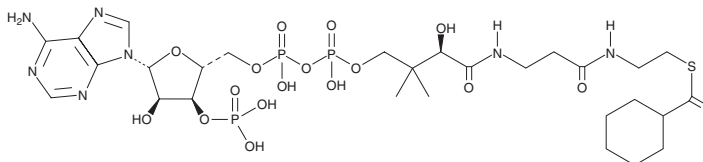
PRODUCT INFORMATION



Cyclohexanoyl-Coenzyme A

Item No. 23504

CAS Registry No.: 5960-12-3
Formal Name: S-cyclohexanecarboxylate
coenzyme A
Synonyms: CHCoA,
Cyclohexanecarboxyl-CoA
MF: $C_{28}H_{46}N_7O_{17}P_3S$
FW: 877.7
Purity: $\geq 95\%$
UV/Vis.: λ_{max} : 260 nm
Supplied as: A crystalline 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

Cyclohexanoyl-Coenzyme A (CHCoA) is supplied as a crystalline solid. Aqueous solutions of CHCoA can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of CHCoA in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

CHCoA is an acyl-CoA that contains a cyclohexane group. It is the activated form of cyclohexane carboxylic acid (CHC) in *R. palustris*.¹ CHC is converted to CHCoA by a succinyl-CoA CHC CoA transferase, and CHCoA is then degraded by a dehydrogenase. CHCoA is converted to hippuric acid in submitochondrial fractions isolated from guinea pig liver.²

References

1. Kung, J.W., Meier, A.-K., Mergelsberg, M., *et al.* Enzymes involved in a novel anaerobic cyclohexane carboxylic acid degradation pathway. *J. Bacteriol.* **196**(20), 3667-3674 (2014).
2. Svardakm A.M. and Scheline, R.R. The aromatization of cyclohexanecarboxyl-CoA to hippuric acid by guinea pig liver mitochondria: Submitochondrial localization. *Mol. Cell. Biochem.* **67**(2), 181-188 (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, 11/15/2022

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