

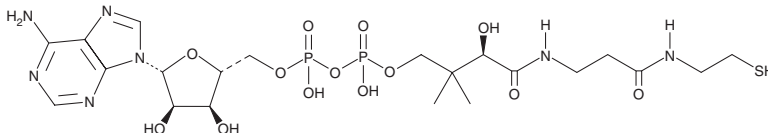
PRODUCT INFORMATION



3'-Dephosphocoenzyme A

Item No. 27390

CAS Registry No.: 3633-59-8
Formal Name: 3'-O-dephosphono-coenzyme A
Synonyms: depCoA, Dephospho-CoA
MF: $C_{21}H_{35}N_7O_{13}P_2S$
FW: 687.6
Purity: $\geq 90\%$
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

3'-Dephosphocoenzyme A is supplied as a crystalline solid. A stock solution may be made by dissolving the 3'-dephosphocoenzyme A in water. The solubility of 3'-dephosphocoenzyme A in water is approximately 50 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

3'-Dephosphocoenzyme A is an intermediate in the biosynthesis of coenzyme A (CoA; Item Nos. 16147 | 21499 | 21722) from pantothenic acid (vitamin B₅; Item No. 17288).¹ It is phosphorylated by CoA synthase in humans to form CoA. 3'-Dephosphocoenzyme A can be used as a transcription initiator in the synthesis of CoA-RNA by *in vitro* transcription.^{2,3}

References

- Leonardi, R., Zhang, Y.M., Rock, C.O., *et al.* Coenzyme A: Back in action. *Prog. Lipid Res.* **44**(2-3), 125-153 (2005).
- Huang, F. Efficient incorporation of CoA, NAD and FAD into RNA by *in vitro* transcription. *Nucleic Acids Res.* **31**(3), e8 (2003).
- Sapkota, K. and Huang, F. Efficient one-pot enzymatic synthesis of dephospho coenzyme A. *Bioorg. Chem.* **76**, 23-27 (2018).

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, 12/05/2022

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