

PRODUCT INFORMATION

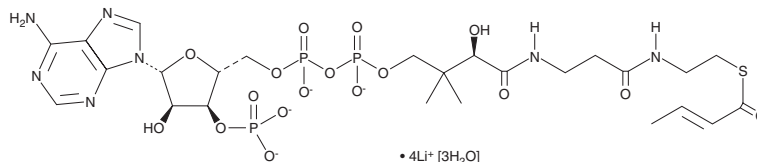


Crotonyl-Coenzyme A (lithium salt hydrate)

Item No. 28606

Formal Name: (2R,3S,4R,5R)-5-(6-amino-9H-purin-9-yl)-2-((((((R)-4-((3-((2-((E)-but-2-enoyl)thio)ethyl)amino)-3-oxopropyl)amino)-3-hydroxy-2,2-dimethyl-4-oxobutoxy)oxidophosphoryl)oxy)oxidophosphoryl)oxy)methyl)-4-hydroxytetrahydrofuran-3-yl phosphate, lithium salt hydrate

Synonym: Crotonyl-CoA
MF: $C_{25}H_{36}N_7O_{17}P_3S \cdot 4Li [3H_2O]$
FW: 913.4
Purity: $\geq 90\%$
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

Crotonyl-coenzyme A (crotonyl-CoA) is supplied as a solid. A stock solution may be made by dissolving the crotonyl-CoA in the solvent of choice, which should be purged with an inert gas. Crotonyl-CoA is slightly soluble (0.1-1 mg/ml) in acetonitrile:water (1:1).

Description

Crotonyl-CoA is a derivative of CoA (Item Nos. 21499 | 21722) that contains the short-chain fatty acid crotonic acid (Item No. 43903).¹ It is formed from lysine and tryptophan catabolism and fatty acid β -oxidation and serves as a precursor to acetyl-CoA. Crotonyl-CoA is a cofactor for crotonyl transferases, which mediate the post-translational modification of lysine residues on various proteins, including histones.² Histone crotonylation is associated with activation of gene expression.

References

1. Wang, Y., Chen, H.-Z., and Tang, X. Crotonyl-coenzyme A (crotonyl-CoA). *Trends in Endocrinol. Metab.* **36**(12), 1154-1155 (2025).
2. Westerveld, M., Besermenji, K., Aidukas, D., et al. Cracking lysine crotonylation (Kcr): Enlightening a promising post-translational modification. *Chembiochem* **26**, e202400639 (2025).

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, 07/01/2026

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