

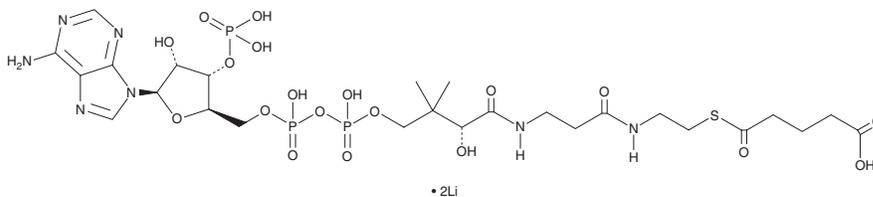
PRODUCT INFORMATION



Glutaryl-Coenzyme A (lithium salt)

Item No. 44592

CAS Registry No.: 103192-48-9
Formal Name: S-(hydrogen pentanedioate) coenzyme A, dilithium salt
Synonym: Glutaryl-CoA
MF: 895.5
FW: $C_{26}H_{42}N_7O_{19}P_3S \cdot 2Li$
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

Glutaryl-coenzyme A (glutaryl-CoA) (lithium salt) is supplied as a solid. A stock solution may be made by dissolving the glutaryl-CoA (lithium salt) in the solvent of choice, which should be purged with an inert gas. Glutaryl-CoA (lithium salt) is sparingly soluble (1-10 mg/ml) in a 1:1 solution of acetonitrile:water.

Description

Glutaryl-CoA is an intermediate in lysine, hydroxylysine, and tryptophan catabolism and a derivative of CoA (Item Nos. 16147 | 21499 | 21722) that contains the short-chain fatty acid glutaric acid.¹ It inhibits α -ketoglutarate dehydrogenase ($K_i = 0.35$ mM).

Reference

1. Sauer, S.W., Okun, J.G., Schwab, M.A., *et al.* Bioenergetics in glutaryl-coenzyme a dehydrogenase deficiency. A role for glutaryl-coenzyme A. *J. Biol. Chem.* **290**(23), 21830-21836 (2005).

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, 02/02/2026

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