

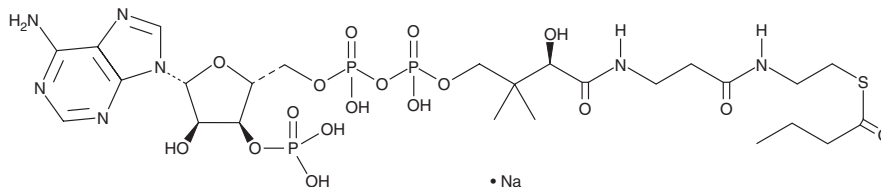
PRODUCT INFORMATION



Butyryl-Coenzyme A (sodium salt)

Item No. 27865

CAS Registry No.: 125527-24-4
Formal Name: S-butanoyl coenzyme A, monosodium salt
Synonyms: S-Butanoyl-CoA, Butyryl-CoA
MF: C₂₅H₄₂N₇O₁₇P₃S • Na
FW: 860.6
Purity: ≥95%
UV/Vis.: λ_{max}: 258 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

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

Butyryl-CoA is a short-chain acyl CoA. It is an intermediate in the synthesis of butyrate, which is produced by colonic bacteria and is important for maintaining the colonic environment.¹

Reference

1. Trachsel, J., Bayles, D.O., Looft, T., *et al.* Function and phylogeny of bacterial butyryl coenzyme A:Acetate transferases and their diversity in the proximal colon of swine. *Appl. Environ. Microbiol.* **82(22)**, 6788-6798 (2016).

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, 10/06/2022

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