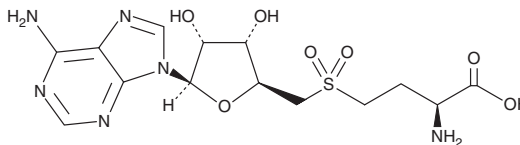


PRODUCT INFORMATION



SAHO₂
Item No. 28726

CAS Registry No.: 53199-56-7
Formal Name: 5'-[[[(3S)-3-amino-3-carboxypromyl]sulfonyl]-5'-deoxy-adenosine
MF: C₁₄H₂₀N₆O₇S
FW: 416.4
Purity: ≥95%
UV/Vis.: λ_{max}: 235 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

SAHO₂ is supplied as a crystalline solid. A stock solution may be made by dissolving the SAHO₂ in the solvent of choice, which should be purged with an inert gas. SAHO₂ is soluble in the organic solvent DMSO at a concentration of approximately 30 mg/ml.

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 SAHO₂ can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of SAHO₂ in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

SAHO₂ is a sulfone form of the methyl donor S-(5'-adenosyl)-L-methionine chloride (SAM; Item No. 13956) and is a substrate for radical SAM enzymes.¹ It is reductively cleaved to 5'-deoxyadenosine (Item No. 29619) and 5'-thioadenosine sulfinic acid by the radical SAM enzymes NosL or NosN.

Reference

1. Mandalapu, D., Ji, X., and Zhang, Q. Reductive cleavage of sulfoxide and sulfone by two radical S-adenosyl-L-methionine enzymes. *Biochemistry* **58(1)**, 36-39 (2019).

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, 03/19/2024

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