

PRODUCT INFORMATION



S-(5'-Adenosyl)-L-methionine (sulfate tosylate)

Item No. 17595

CAS Registry No.: 97540-22-2

Formal Name: 5'-[[[(3S)-3-amino-3-carboxypropyl]methylsulfonio]-5'-deoxy-adenosine, monosulfate, 4-methylbenzenesulfonate monosulfate

Synonyms: AdoMet, FO 1561, SAM, SAME

MF: $C_{15}H_{23}N_6O_5S \cdot HSO_4 \cdot H_2SO_4 \cdot C_7H_8O_3S$

FW: 766.8

Purity: $\geq 95\%$

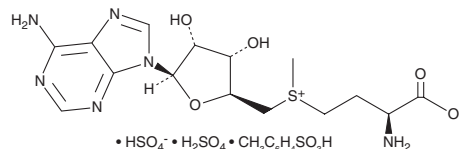
UV/Vis.: λ_{max} : 260 nm

Supplied as: A solid

Storage: $-20^\circ C$

Stability: ≥ 4 years

Special Conditions: Hygroscopic



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

S-(5'-Adenosyl)-L-methionine (SAM) (sulfate tosylate) is supplied as a solid. A stock solution may be made by dissolving the SAM (sulfate tosylate) in the solvent of choice, which should be purged with an inert gas. SAM (sulfate tosylate) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of SAM (sulfate tosylate) in ethanol is approximately 10 mg/ml and approximately 30 mg/ml in DMSO and DMF.

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

Description

SAM is a methyl donor and cofactor.¹ It is involved in several biological processes, including epigenetic regulation, post-translational modification of proteins, and metabolism. Formulations containing SAM have been used as dietary supplements.

Reference

1. Loenen, W.A.M. S-Adenosylmethionine: Jack of all trades and master of everything? *Biochem. Soc. Trans.* **34**(2), 330-333 (2006).

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, 09/06/2023

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