

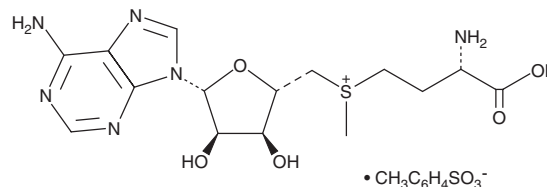
PRODUCT INFORMATION



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

Item No. 16376

CAS Registry No.: 52248-03-0
Formal Name: 5'-[[[(3S)-3-amino-3-carboxypropyl]methylsulfonio]-5'-deoxy-adenosine, 4-methylbenzenesulfonate
Synonym: AdoMet, SAM, SAMe
MF: C₁₅H₂₃N₆O₅S • C₇H₇O₃S
FW: 570.6
Purity: ≥90%
UV/Vis.: λ_{max}: 221, 262 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

S-(5'-Adenosyl)-L-methionine (SAM) (tosylate) is supplied as a crystalline solid. A stock solution may be made by dissolving the SAM (tosylate) in the solvent of choice, which should be purged with an inert gas. SAM (tosylate) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of SAM (tosylate) in these solvents is approximately 10 and 5 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 SAM (tosylate) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of SAM (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, 10/26/2022

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