

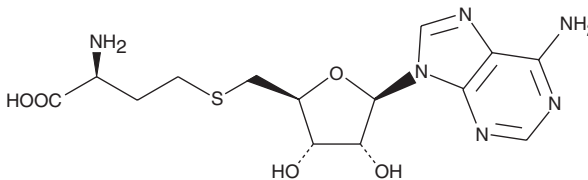
PRODUCT INFORMATION



S-Adenosylhomocysteine

Item No. 13603

CAS Registry No.: 979-92-0
Formal Name: S-(5'-deoxyadenosin-5'-yl)-L-homocysteine
Synonyms: AdoHcy, SAH
MF: C₁₄H₂₀N₆O₅S
FW: 384.4
Purity: ≥98%
UV/Vis.: λ_{max}: 260 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-Adenosylhomocysteine is supplied as a crystalline solid. A stock solution may be made by dissolving the S-adenosylhomocysteine in the solvent of choice, which should be purged with an inert gas. S-Adenosylhomocysteine is soluble in the solvent 1M HCl at a concentration of approximately 20 mg/ml.

Description

S-Adenosyl-L-homocysteine (SAH) is an amino acid derivative and an intermediate, by-product, or modulator of several metabolic pathways, including the activated methyl cycle and cysteine biosynthesis. It is also a product of S-adenosyl-methionine (SAM)-dependent methylation of biological molecules, including DNA, RNA, and histones and other proteins. SAH is a risk factor for many diseases, including cancer and neurodegenerative diseases. In addition, inhibitors that block its hydrolysis are being developed as anti-viral, anti-parasitic, anti-arthritis and immunosuppressive agents.

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, 11/18/2022

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