

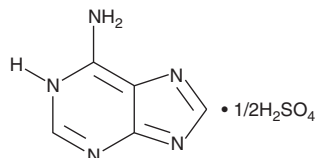
PRODUCT INFORMATION



Adenine (hemisulfate)

Item No. 18148

CAS Registry No.: 321-30-2
Formal Name: 9H-purine-6-amine, hemisulfate
MF: C₅H₅N₅ • 1/2H₂O₄S
FW: 184.2
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

Adenine (hemisulfate) is supplied as a crystalline solid. A stock solution may be made by dissolving the adenine (hemisulfate) in the solvent of choice, which should be purged with an inert gas. Adenine (hemisulfate) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of adenine (hemisulfate) in these solvents is approximately 30 and 2 mg/ml, respectively.

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 adenine (hemisulfate) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of adenine (hemisulfate) in PBS, pH 7.2, is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Adenine is a purine base that acts as a precursor, substrate, or cofactor in diverse biochemical pathways.¹ It is a fundamental component of DNA, RNA, and ATP, as well as numerous other natural biochemicals.

Reference

1. Nuki, G. Human purine metabolism: Some recent advances and relationships with immunodeficiency. *Ann. Rheum. Dis.* **42**, 8-11 (1983).

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, 12/22/2022

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