

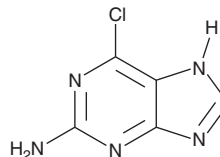
PRODUCT INFORMATION



2-Amino-6-chloropurine

Item No. 30400

CAS Registry No.: 10310-21-1
Formal Name: 6-chloro-9H-purin-2-amine
Synonyms: 6-Chloroguanine, NSC 29570
MF: C₅H₄ClN₅
FW: 169.6
Purity: ≥98%
UV/Vis.: λ_{max}: 218 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

2-Amino-6-chloropurine is supplied as a solid. A stock solution may be made by dissolving the 2-amino-6-chloropurine in the solvent of choice, which should be purged with an inert gas. 2-Amino-6-chloropurine is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of 2-amino-6-chloropurine in these solvents is approximately 3 mg/ml.

2-Amino-6-chloropurine is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 2-amino-6-chloropurine should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. 2-Amino-6-chloropurine has a solubility of approximately 0.12 mg/ml in a 1:7 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

2-Amino-6-chloropurine is a precursor in the synthesis of nucleoside analogs with antiviral activity against Epstein-Barr virus (EBV) and human herpes virus 6 (HHV-6).¹

Reference

1. Qiu, Y.-L., Ksebati, M.B., Ptak, R.G., *et al.* (Z)- and (E)-2-((hydroxymethyl)cyclopropylidene)methyladenine and -guanine. New nucleoside analogues with a broad-spectrum antiviral activity. *J. Med. Chem.* **41(1)**, 10-23 (1998).

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, 05/18/2020

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