

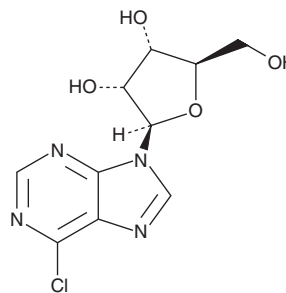
PRODUCT INFORMATION



6-Chloropurine Riboside

Item No. 36002

CAS Registry No.: 5399-87-1
Formal Name: 6-chloro-9-β-D-ribofuranosyl-9H-purine
Synonym: NSC 4910
MF: C₁₀H₁₁ClN₄O₄
FW: 286.7
Purity: ≥98%
Supplied as: A 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

6-Chloropurine riboside is supplied as a solid. A stock solution may be made by dissolving the 6-chloropurine riboside in the solvent of choice, which should be purged with an inert gas. 6-Chloropurine riboside is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of 6-chloropurine riboside in these solvents is approximately 5 and 2 mg/ml, respectively.

Description

6-Chloropurine riboside is a nucleoside precursor.¹ It has been used in the synthesis of adenosine derivatives with antiproliferative activity in human gastric cancer cells expressing the adenosine A₃ receptor. 6-Chloropurine riboside has also been used in the synthesis of N⁶-benzyladenosine derivatives with antiproliferative activity in a variety of cancer cells or inhibitory activity against *T. gondii* replication in CRL-1634/Hs27 human skin fibroblasts.^{2,3}

References

1. Zurita, F.V., Vega, N.B., and Cabrera, M.G. Semisynthesis, characterization and evaluation of new adenosine derivatives as antiproliferative agents. *Molecules* **23**(5), 1111 (2018).
2. Dolezal, K., Popa, I., Hauserová, E., *et al.* Preparation, biological activity and endogenous occurrence of N⁶-benzyladenosines. *Bioorg. Med. Chem.* **15**(11), 3737-3747 (2007).
3. Kim, Y.A., Sharon, A., Chu, C.K., *et al.* Synthesis, biological evaluation and molecular modeling studies of N⁶-benzyladenosine analogues as potential anti-toxoplasma agents. *Biochem. Pharmacol.* **73**(10), 1558-1572 (2007).

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, 12/09/2022

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