

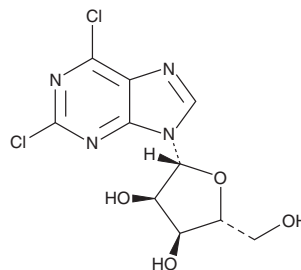
PRODUCT INFORMATION



2,6-Dichloropurine-9-β-D-ribose

Item No. 31405

CAS Registry No.: 13276-52-3
Formal Name: 2,6-dichloro-9-β-D-ribofuranosyl-9H-purine
Synonym: 2,6-Dichloropurine riboside
MF: C₁₀H₁₀Cl₂N₄O₄
FW: 321.1
Purity: ≥98%
UV/Vis.: λ_{max}: 274 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

2,6-Dichloropurine-9-β-D-ribose is supplied as a crystalline solid. A stock solution may be made by dissolving the 2,6-dichloropurine-9-β-D-ribose in the solvent of choice, which should be purged with an inert gas. 2,6-Dichloropurine-9-β-D-ribose is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of 2,6-dichloropurine-9-β-D-ribose in these solvents is approximately 30 mg/ml.

2,6-Dichloropurine-9-β-D-ribose is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 2,6-dichloropurine-9-β-D-ribose should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. 2,6-Dichloropurine-9-β-D-ribose has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

2,6-Dichloropurine-9-β-D-ribose is a building block.^{1,2} It has been used in the synthesis of photoaffinity probes for nucleotide binding sites in proteins.

References

1. Wower, J., Hixson, S.S., Sylvers, L.A., *et al.* Synthesis of 2,6-diazido-9-(β-D-ribofuranosyl)purine 3',5'-bisphosphate: Incorporation into transfer RNA and photochemical labeling of *Escherichia coli* ribosomes. *Bioconjug. Chem.* **5**(2), 158-161 (1994).
2. Seiter, M.S., Bauer, M.P., Bogel, P.D., *et al.* Synthesis of novel spin-labeled photoaffinity derivatives of NAD⁺ and ATP and their characterization as coenzymes and substrates of several enzymes. *Synthesis* **2**, 269-273 (1996).

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