

PRODUCT INFORMATION



6-Chloropurine riboside-5'-O-diphosphate (sodium salt)

Item No. 38552

Formal Name: 6-chloro-9-[5-O-[hydroxy(phosphonooxy)phosphinyl]-β-D-ribofuranosyl]-9H-purine, trisodium salt

Synonym: 6-chloro PuTP

MF: C₁₀H₁₀ClN₄O₁₀P₂ • 3Na

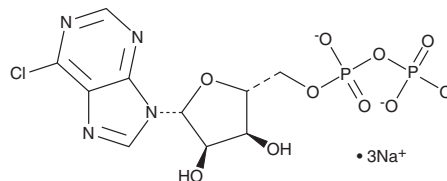
FW: 512.6

Purity: ≥95%

Supplied as: A solution in water

Storage: -80°C

Stability: ≥4 years



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

Description

6-Chloropurine riboside-5'-O-diphosphate is a synthetic precursor and phosphorylated form of 6-chloropurine riboside (Item No. 36002).^{1,2} It has been used in the synthesis of cytokinins with anticancer activity and spin-labeled derivatives of NAD⁺ (Item No. 16077).

References

1. Voller, J., Zatloukal, M., Lenobel, R., *et al.* Anticancer activity of natural cytokinins: A structure-activity relationship study. *Phytochemistry* **71**(11-12), 1350-1359 (2010).
2. Wenzel, H.R., Pfeleiderer, G., Trommer, W.E., *et al.* The synthesis of spin-label derivatives of NAD⁺ and its structural components and their binding to lactate dehydrogenase. *Biochim. Biophys. Acta* **452**(2), 292-301 (1976).

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, 04/23/2026

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