

PRODUCT INFORMATION



Isoguanosine-5'-O-diphosphate (sodium salt)

Item No. 38566

Formal Name: ((2R,3S,4R,5R)-5-(6-amino-2-oxo-2,3-dihydro-9H-purin-9-yl)-3,4-dihydroxytetrahydrofuran-2-yl)methyl diphosphate, trisodium salt
Synonym: isoGDP

MF: $C_{10}H_{12}N_5O_{11}P_2 \cdot 3Na$

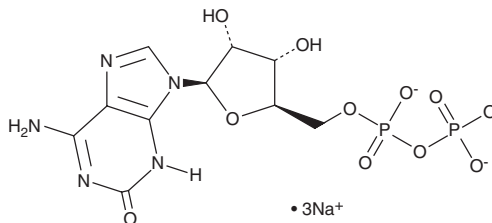
FW: 509.1

Purity: $\geq 95\%$

Supplied as: A solution in water

Storage: $-80^{\circ}C$

Stability: ≥ 4 years



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

Description

Isoguanosine-5'-O-diphosphate (isoGDP) is an isomer of GDP (Item No. 31704) and a phosphorylated form of crotonoside (Item No. 28219). isoGDP is an inhibitor of mitochondrial ATPase ($IC_{50} = 3 \mu M$).¹

Reference

1. Harris, D.A., Gomez-Fernandez, J.C., Klungsøyr, L., *et al.* Specificity of nucleotide binding and coupled reactions utilising the mitochondrial ATPase. *Biochim. Biophys. Acta* **504(3)**, 364-383 (1978).

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