

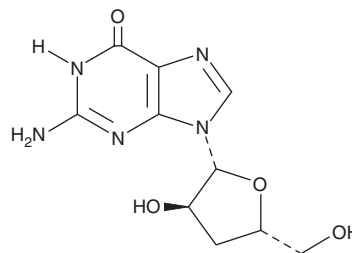
# PRODUCT INFORMATION



## 3'-Deoxyguanosine

Item No. 17907

CAS Registry No.: 3608-58-0  
Formal Name: 3'-deoxy-guanosine  
MF:  $C_{10}H_{13}N_5O_4$   
FW: 267.2  
Purity:  $\geq 95\%$   
UV/Vis.:  $\lambda_{\max}$ : 254 nm  
Supplied as: A crystalline solid  
Storage:  $-20^{\circ}\text{C}$   
Stability:  $\geq 4$  years



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

### Laboratory Procedures

3'-Deoxyguanosine is supplied as a crystalline solid. A stock solution may be made by dissolving the 3'-deoxyguanosine in the solvent of choice, which should be purged with an inert gas. 3'-Deoxyguanosine is soluble in the organic solvent DMSO at a concentration of approximately 10 mg/ml. The solubility of 3'-deoxyguanosine in 0.1 M HCl is approximately 10 mg/ml.

### Description

3'-Deoxyguanosine is a ligand that can be complexed with enzymes, such as purine nucleoside phosphorylase, and receptors, in order to study structure-activity relationships.<sup>1,2</sup> This compound can also be used to selectively impair transcription.<sup>3</sup>

### References

1. Canduri, F., Silva, R.G., dos Santos, D.M., *et al.* Structure of human PNP complexed with ligands. *Acta Crystallogr. D Biol. Crystallogr.* **61**(Pt 7), 856-862 (2005).
2. Traversa, U., Bombi, G., Camaioni, E., *et al.* Rat brain guanosine binding site. Biological studies and pseudo-receptor construction. *Bioorg. Med. Chem.* **11**(24), 5417-5425 (2003).
3. King, R.W., Zecher, M., and Jefferies, M.W. Inhibition of the replication of a hepatitis C virus-like RNA template by interferon and 3'-deoxycytidine. *Antivir. Chem. Chemother.* **13**(6), 363-370 (2002).

#### 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, 11/10/2022

#### 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