

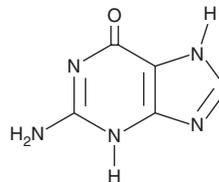
PRODUCT INFORMATION



Guanine

Item No. 34248

CAS Registry No.: 73-40-5
Formal Name: 2-amino-1,9-dihydro-6H-purin-6-one
MF: C₅H₅N₅O
FW: 151.1
Purity: ≥95%
UV/Vis.: λ_{max}: 248 nm
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

Guanine is supplied as a solid. A stock solution may be made by dissolving the guanine in water or aqueous acid. Guanine is soluble in aqueous acid and sparingly soluble in water. We do not recommend storing the aqueous solution for more than one day.

Description

Guanine is a purine base.¹ It forms complementary base pairs with the pyrimidine cytosine in DNA and RNA. Guanine can be linked to ribose via a β-N₉-glycosidic bond to form the purine nucleoside guanosine (Item No. 27702), which can be phosphorylated to guanine-based purines that have numerous roles in intracellular signal transduction.^{2,3}

References

1. Berg, J.M., Tymoczko, J.L., and Stryer, L. *Biochemistry*. 5th ed., W.H. Freeman, New York (2002).
2. Devereaux, Z.J., He, C.C., Zhu, Y., *et al.* Structures and relative glycosidic bond stabilities of protonated 2'-fluoro-substituted purine nucleosides. *J. Am. Soc. Mass Spectrom.* **30(8)**, 1521-1536 (2019).
3. Di Liberto, V., Mudò, G., Garozzo, R., *et al.* The guanine-based purinergic system: The tale of an orphan neuromodulation. *Front. Pharmacol.* **7**, 158 (2016).

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, 11/08/2022

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