

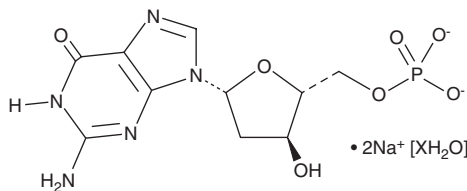
PRODUCT INFORMATION



2'-Deoxyguanosine 5'-monophosphate (sodium salt hydrate)

Item No. 16408

CAS Registry No.: 146877-98-7
Formal Name: 2'-deoxy-5'-guanylic acid, disodium salt hydrate
Synonym: dGMP
MF: C₁₀H₁₂N₅O₇P • 2Na [XH₂O]
FW: 391.2
Purity: ≥95%
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'-Deoxyguanosine 5'-monophosphate (dGMP) (sodium salt hydrate) is supplied as a crystalline solid. Aqueous solutions of dGMP (sodium salt hydrate) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of dGMP (sodium salt hydrate) in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

dGMP is used as a substrate of guanylate kinases to generate dGDP, which in turn is phosphorylated to dGTP, a nucleotide precursor used in DNA synthesis. In addition, dGMP can be efficiently phosphorylated by the type I thymidylate kinase of the malarial agent *P. falciparum*.^{1,2} In humans, deficiency of mitochondrial deoxyguanosine kinase, which synthesizes dGMP from ATP and deoxyguanosine, contributes to mitochondrial DNA depletion syndrome.^{3,4}

References

1. Kandeel, M., Kitamura, Y., and Kitade, Y. The exceptional properties of Plasmodium deoxyguanylate pathways as a potential area for metabolic and drug discovery studies. *Nucleic Acids Symp. Ser. (Oxf)* **53**, 39-40 (2009).
2. Whittingham, J.L., Carrero-Lerida, J., Brannigan, J.A., et al. Structural basis for the efficient phosphorylation of AZT-MP (3'-azido-3'-deoxythymidine monophosphate) and dGMP by Plasmodium falciparum type I thymidylate kinase. *Biochem. J.* **428(3)**, 499-509 (2014).
3. Taanman, J.W., Muddle, J.R., and Muntau, A.C. Mitochondrial DNA depletion can be prevented by dGMP and dAMP supplementation in a resting culture of deoxyguanosine kinase-deficient fibroblasts. *Hum. Mol. Genet.* **12(15)**, 1839-1845 (2003).
4. Bulst, S., Abicht, A., Holinski-Feder, E., et al. In vitro supplementation with dAMP/dGMP leads to partial restoration of mtDNA levels in mitochondrial depletion syndromes. *Hum. Mol. Genet.* **18(9)**, 1590-1599 (2009).

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, 12/15/2022

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