

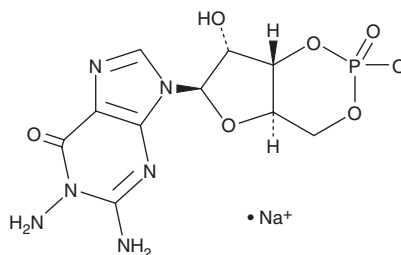
PRODUCT INFORMATION



1-Aminoguanosine-3',5'-monophosphate (sodium salt)

Item No. 38638

CAS Registry No.: 1380911-23-8
Formal Name: 1-amino-cyclic 3',5'-(hydrogen phosphate) guanosine, monosodium salt
Synonym: 1-amino cGMP
MF: C₁₀H₁₂N₆O₇P • Na
FW: 382.2
Purity: ≥95%
Supplied as: A solid
Storage: -80°C
Stability: ≥2 years



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

Laboratory Procedures

1-Aminoguanosine-3',5'-monophosphate (1-amino cGMP) (sodium salt) is supplied as a solid. A stock solution may be made by dissolving the 1-amino cGMP (sodium salt) in water. We do not recommend storing the aqueous solution for more than one day.

Description

1-amino cGMP is a derivative of the secondary messenger cGMP (Item No. 18821) and a substrate for cGMP-dependent protein kinase.¹

Reference

1. Corbin, J.D., Og Reid, D., Miller, J.P., *et al.* Studies of cGMP analog specificity and function of the two intrasubunit binding sites of cGMP-dependent protein kinase. *J. Biol. Chem.* **261(3)**, 1208-1214 (1986).

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, 08/08/2023

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