

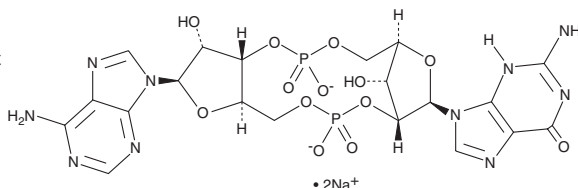
PRODUCT INFORMATION



2'3'-cGAMP (sodium salt)

Item No. 19887

CAS Registry No.: 2734858-36-5
Formal Name: adenylyl-(3'→5')-2'-guanylic acid, cyclic nucleotide, disodium salt
Synonyms: 2'3'-Cyclic GMP-AMP, Guanosine-Adenosine 2',3'-cyclic monophosphate
MF: C₂₀H₂₂N₁₀O₁₃P₂ • 2Na
FW: 718.4
Purity: ≥98%
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

2'3'-cGAMP (sodium salt) is supplied as a solid. A stock solution may be made by dissolving the 2'3'-cGAMP (sodium salt) in water. We do not recommend storing the aqueous solution for more than one day.

Description

2'3'-cGAMP is a second messenger produced from ATP and GTP by cGMP-AMP synthase (cGAS) in the cytoplasm of mammalian cells in response to the presence of DNA.¹⁻³ It contains a non-canonical 2'5'-phosphodiester bond and binds tightly to the adapter protein STING (stimulator of interferon genes) with a K_d value of 3.79 nM, which is significantly lower than that observed for 2'2'-cGAMP, 3'3'-cGAMP (Item No. 17966), cyclic di-GMP (Item No. 17144), and 3'2'-cGAMP (K_ds = 0.287, 1.04, 1.21, and 1.61 μM, respectively).³ It is a potent inducer of IFN-β mRNA expression with an EC₅₀ value of 19.4 nM in L929 cells.³

References

1. Wu, J., Sun, L., Chen, X., *et al.* Cyclic GMP-AMP is an endogenous second messenger in innate immune signaling by cytosolic DNA. *Science* **339**(6121), 826-830 (2013).
2. Sun, L., Wu, J., Du, F., *et al.* Cyclic GMP-AMP synthase is a cytosolic DNA sensor that activates the type I interferon pathway. *Science* **339**(6121), 786-791 (2013).
3. Zhang, X., Shi, H., Wu, J., *et al.* Cyclic GMP-AMP containing mixed phosphodiester linkages is an endogenous high-affinity ligand for STING. *Mol. Cell* **51**(2), 226-235 (2015).

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, 02/10/2025

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