

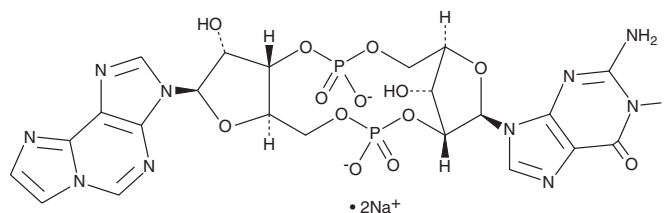
PRODUCT INFORMATION



ϵ -2'3'-cGAMP (sodium salt)

Item No. 28321

Formal Name: N,1-(1,2-ethenediyl)adenylyl-(3'→5')-2'-guanylic acid, cyclic nucleotide, disodium salt
MF: $C_{22}H_{22}N_{10}O_{13}P_2 \cdot 2Na$
FW: 742.4
Purity: $\geq 95\%$
Ex./Em. Max: 300/408 nm
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



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

Laboratory Procedures

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

Description

ϵ -2'3'-cGAMP is a fluorescent derivative of the second messenger 2'3'-cGAMP (Item No. 19887). It has activity in an interferon-stimulated genes (ISG) reporter assay using THP-1 cells ($EC_{50} = \leq 100$ nM).¹ ϵ -2'3'-cGAMP displays excitation/emission maxima of 300/408 nm, respectively.

Reference

1. Zhou, B., Sun, L., Wei, Q., *et al.* Cyclic di-nucleotide compounds and methods of use. *Immune Sensor, LLC*. WO2017/161349A1 (2017).

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, 09/14/2023

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