

# pppApG

## Potential endogenous metabolite of the metazoan second messenger 2',3'-cGAMP and/or of the bacterial signalling nucleotide 3',3'-cGAMP

In pppApG a 5'-ATP unit is connected with a 5'-GMP unit via a 3' → 5' linkage to form a linear dinucleotide. pppApG is a potential endogenous metabolite of the metazoan second messenger 2',3'-cGAMP (c[G(2',5')pA(3',5')p], Prod. No. BLG-C161) and/or of the bacterial signalling nucleotide 3',3'-cGAMP (c-(ApGp), Prod. No. BLG-C117). pppApG was reported to be the linear dinucleotide intermediate in the enzymatic production of 3',3'-cGAMP by the *Vibrio cholerae* pathogenicity factor DncV.

*BLG-T050-05 (5 x 1 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.*

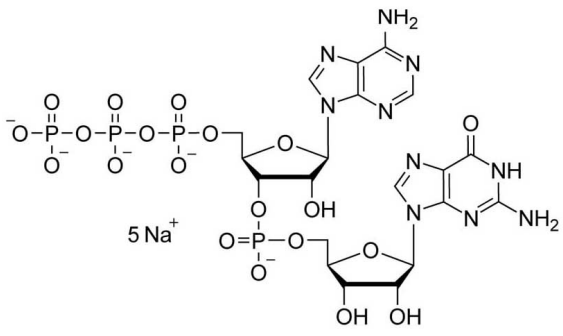
### Ordering Information

[Order Online »](#)

BLG-T050-01	1µmol
BLG-T050-05	5x1µmol

### Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** 5'-Triphosphoadenylyl-(3'→5')-guanosine . sodium salt

**Appearance** 10 mM aqueous solution

**CAS** 68892-47-7

**Couple Type** Modified nucleotides

**Excitation Maximum** 256 nm

**Formula**  $C_{20}H_{28}N_{10}O_{20}P_4$  (free acid)

**MW** 852.39 (free acid)

**Purity** > 95% HPLC

**Technical Info / Product Notes** For the Original Manufacturer's data sheet please [click here](#).



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)