

# Di-MANT-2'3'-cGAMP

## Fluorescent analogue of the metazoan second messenger 2'3'-cGAMP

Potentially membrane permeant, fluorescent analogue of the metazoan cyclic dinucleotide second messenger and STING activator c[G(2',5')pA(3',5')p] (aka cGAMP(2'-5') or 2'3'-cGAMP, Cat. No. BLG-C161), where both, the 2'- and the 3'-hydroxyl group at the ribose moieties, are esterified by the fluorescent MANT group. The fluorophore ( $\lambda_{exc}$  355 nm;  $\lambda_{em}$  448 nm) has a certain sensitivity for its environment and can change its spectral properties upon binding.

*BLG-C284-005 (5 x 0.1  $\mu$ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.*

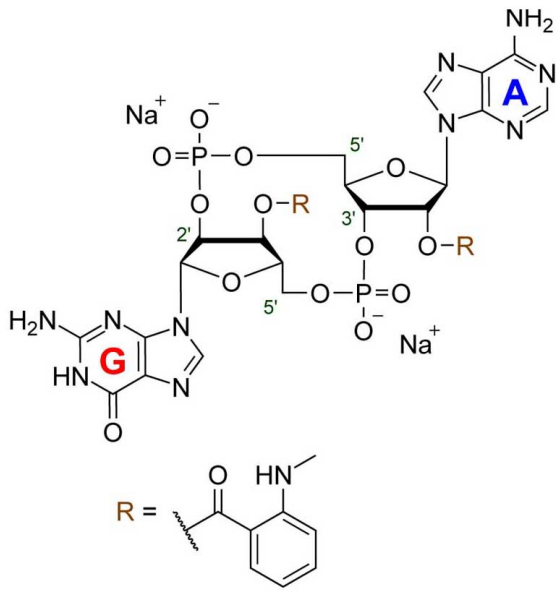
### Ordering Information

[Order Online »](#)

BLG-C284-001	0.1 $\mu$ mol
BLG-C284-005	5x0.1 $\mu$ mol

### Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** c[3'-MANT-G(2',5')p-2'-MANT-A(3',5')p], Cyclic (3'-O-(N'-methylantraniloyl)guanosine-(2'→5')-monophosphate-2'-O-(N'-methylantraniloyl)adenosine-(3'→5')-monophosphate) . sodium salt

**Appearance** Solid

**Formula**  $C_{36}H_{38}N_{12}O_{15}P_2$  (free acid)

**MW** 940.7 (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)