

# c-triGMP

c-triGMP is a cyclic nucleotide in which three 5'-GMP units are interconnected via 3'-5' phosphodiester bonds to form a cyclic structure. c-triGMP is structurally related to c-triAMP (Cat. No. C 362) which is involved in Type III CRISPR-Cas systems in prokaryotes. Reference: Ross et al., J. Biol. Chem., 265, 18933 - 18943 (1990).

*BLG-C367-025 (5 x 0.5 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.*

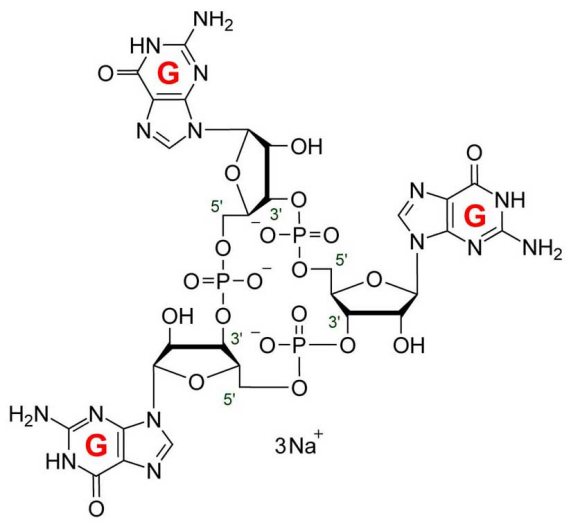
## Ordering Information

[Order Online »](#)

BLG-C367-005	0.5µmol
BLG-C367-025	5x0.5µmol

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Long Term Storage -20°C

Shipping Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** Cyclic triguanosine monophosphate . sodium salt, c[G(3',5')pG(3',5')pG(3',5')p], cyclic triguanylate, c-G3

**CAS** 73353-13-6

**Couple Type** Modified nucleotides

**Formula**  $C_{30}H_{36}N_{15}O_{21}P_3$  (free acid)

**MW** 1035.6 (free acid)

**Purity** > 95% HPLC

**Quantity** 0.5µmol is equivalent to ~0.5 mg

**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)