

# Mant-cGMP

## Fluorescent, PKG-inactive cyclic GMP analogue

Fluorescent, membrane permeant cyclic GMP analogue, suitable for phosphodiesterase studies. The compound shows a decrease in fluorescence when cleaved by PDE.

*BLG-M009-50 (5 x 10  $\mu$ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.*

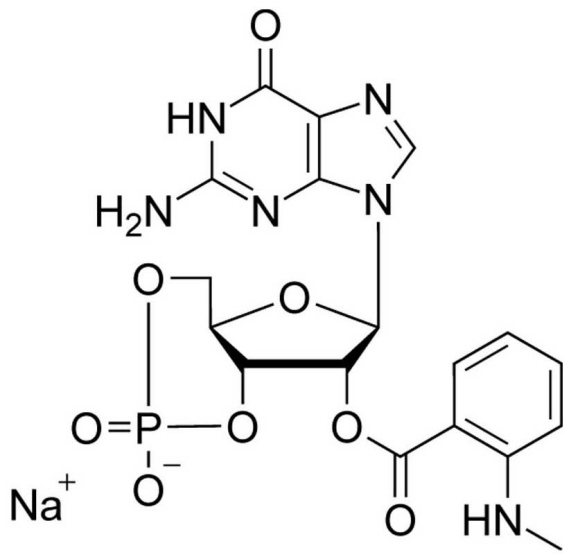
### Ordering Information

[Order Online »](#)

BLG-M009-10	10 $\mu$ mol
BLG-M009-50	5x10 $\mu$ 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** 2'-O-(N<sup>1</sup>-Methylantraniloyl)guanosine-3',5'-cyclic monophosphate . sodium salt

**CAS** 83707-15-7

**Couple Type** Modified nucleotides

**Formula** C<sub>18</sub>H<sub>18</sub>N<sub>6</sub>O<sub>8</sub>P . Na

**MW** 500.3

**Purity** > 98% HPLC

**Quantity** 10µmol≈5mg

**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-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

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

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

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