

# cGMP

## Second messenger

Target receptors for the second messenger cyclic GMP include protein kinase G isozymes, cGMP-dependent ion channels as well as cGMP-responsive phosphodiesterases. Cyclic GMP is metabolized rapidly in presence of corresponding phosphodiesterases.

*BLG-G001-500 (5 x 100  $\mu$ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.*

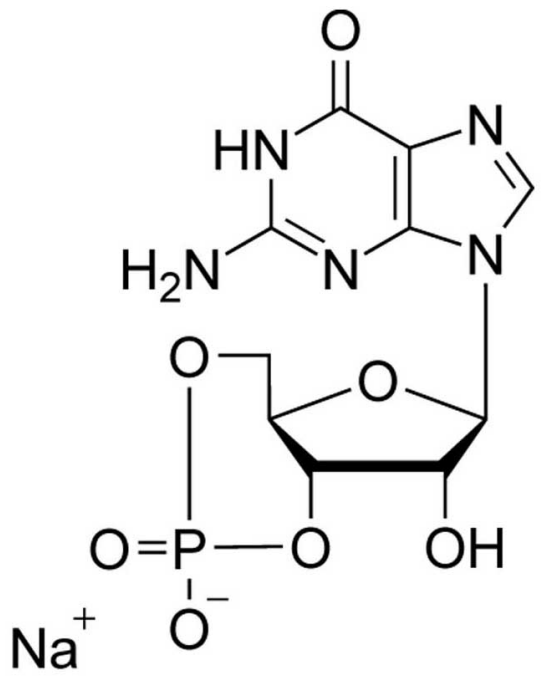
### Ordering Information

[Order Online »](#)

BLG-G001-500	5x100 $\mu$ mol
--------------	-----------------

### 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** Guanosine-3',5'-cyclic monophosphate . sodium salt

**Appearance** Crystallized or lyophilized sodium salt.

**CAS** 40732-48-7

**Couple Type** Activator, Modified nucleotides

**Formula**  $C_{10}H_{11}N_5O_7P \cdot Na$

**MW** 367.2

**Purity** > 98% HPLC

**Quantity** 100µmol≈37mg

**Solubility** Soluble in water or buffer.

**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)