

# Rp-cGMPS

## PDE-resistant PKG inhibitor

Competitive inhibitor of protein kinase G type I alpha (cyclic GMP antagonist). Resistant against mammalian cyclic nucleotide- dependent phosphodiesterases, no metabolic side effects. Both membrane permeability and kinase specificity of this analog are low. Rp-8-pCPT-cGMPS (Prod. No. [BLG-C013](#)) and Rp-8-Br-PET-cGMPS (Prod. No. [BLG-P007](#)) are much more permeant and selective antagonists!

*BLG-G016-25 (5 x 5  $\mu$ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.*

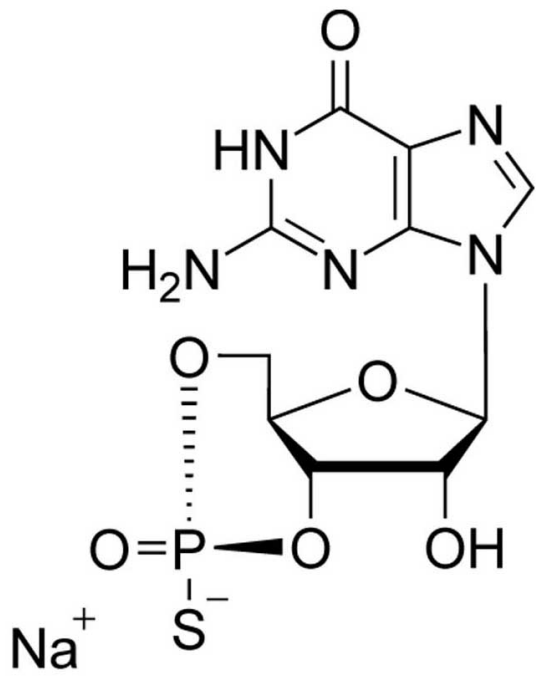
### Ordering Information

[Order Online »](#)

<b>BLG-G016-05</b>	5 $\mu$ mol
<b>BLG-G016-25</b>	5x5 $\mu$ 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 monophosphorothioate . sodium salt, Rp-isomer

**Appearance** Lyophilized or crystallized sodium salt.

**CAS** 86562-09-6

**Couple Type** Modified nucleotides

**Formula**  $C_{10}H_{11}N_5O_6PS \cdot Na$

**MW** 383.3

**Purity** > 99% HPLC

**Quantity** 5µmol≈1.9mg

**Reconstitution** Rinse tube walls carefully and preferably use ultrasonic or vortex to achieve total and uniform mixing.

**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)