

# Rp-8-Br-cGMPS

PDE-resistant inhibitor of protein kinase G I  
 $\alpha$  and I  $\beta$

Competitive inhibitor of protein kinase G type I alpha and I beta (cyclic GMP antagonist). Resistant against mammalian cyclic nucleotide-dependent phosphodiesterases, no metabolic side effects. Significantly more lipophilic compared to cyclic GMP or Rp-cGMPS.

*BLG-B005-25 (5 x 5  $\mu$ mol pack size) is not sold in the U.S. or Canada.  
Please [contact us](#) for available options.*

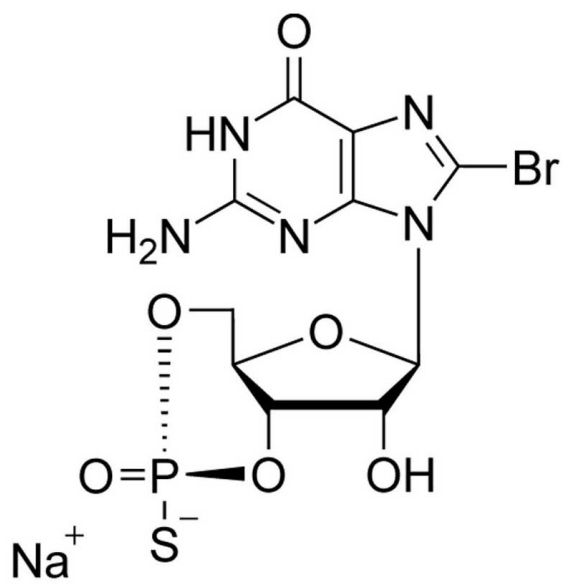
## Ordering Information

[Order Online »](#)

BLG-B005-05	5 $\mu$ mol
BLG-B005-25	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** 8-Bromoguanosine-3',5'-cyclic monophosphorothioate . sodium salt, Rp-isomer

**Appearance** Lyophilized or crystallized sodium salt.

**CAS** 150418-07-8

**Couple Type** Modified nucleotides

**Formula**  $C_{10}H_{10}BrN_5O_6PS \cdot Na$

**MW** 462.2

**Purity** > 99% HPLC

**Quantity** 5µmol≈2.3mg

**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** For the Original Manufacturer's data sheet please [click here](#).

### Notes



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)