

# Rp-8-pCPT-cGMPS

## Membrane-permeant and PDE-resistant PKG inhibitor

Potent selective inhibitor of protein kinase G I  $\alpha$ , I  $\beta$  and especially of type II. Activates retinal cGMP channels and thus discriminates between kinase and channel effects. Absolutely resistant against mammalian cyclic nucleotide-dependent phosphodiesterases, no metabolic side effects. Significantly more lipophilic and membrane permeant compared to Rp-cGMPS (Prod. No. [BLG-G016](#)) or Rp-8-Br-cGMPS (Prod. No. [BLG-B005](#)).

*BLG-C013-05 (5 x 1  $\mu$ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.*

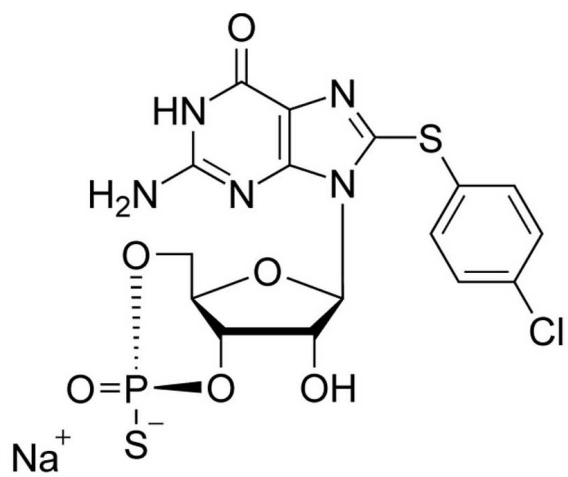
### Ordering Information

[Order Online »](#)

BLG-C013-01	1 $\mu$ mol
BLG-C013-05	5x1 $\mu$ 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	8-(4-Chlorophenylthio)guanosine 3',5'-cyclic monophosphorothioate . sodium salt, Rp-isomer
Appearance	Lyophilized or crystallized sodium salt.
CAS	153660-04-9
Couple Type	Inhibitor, Modified nucleotides
Formula	$C_{16}H_{14}ClN_5O_6PS_2 \cdot Na$
MW	525.9
Purity	> 99% HPLC
Quantity	1µmol≈0.5mg
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 <a href="#">click here</a> .



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)