

# Rp-8-Br-cGMPS

## PKG inhibitor

Cell-permeable, phosphodiesterase-resistant cGMP antagonist. Reversed NO stimulation of random migration in rabbit neutrophils. Used at 30µM to competitively inhibit 8-Br-cGMP relaxation of rabbit aortas. The product is not sterile. Resistant to nucleotide-dependent phosphodiesterases; no metabolic side effects. Significantly more lipophilic and membrane permeant as compared to cyclic GMP or Rp-cGMPS.

Citations: 5

[View Online »](#)

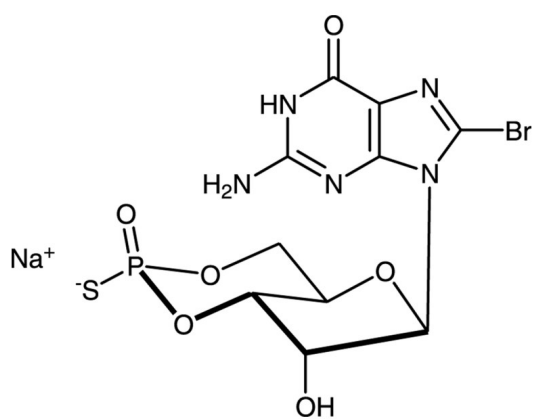
## Ordering Information

[Order Online »](#)

BML-CN216-0500	500µg
----------------	-------

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, at -20°C for up to 1 year. Store solutions at -20°C for up to 3 months.
Long Term Storage	-20°C
Shipping	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

Appearance	White lyophilized powder.
CAS	150418-07-8
Couple Target	PKG
Couple Type	Inhibitor
Formula	$C_{10}H_{10}BrN_5O_6PS \cdot Na$
MW	462.2
Purity	≥98% (HPLC)
Solubility	Soluble in water (25mg/ml).

Last modified: May 29, 2024



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)