

# 8-Br-PET-cGMP

## Site-selective analogue of cyclic GMP

Potent membrane permeant activator of cGMP-dependent protein kinase type I  $\alpha$  and I  $\beta$ . In contrast to common stimulators such as 8-Br-cGMP or 8-pCPT-cGMP which mainly activate kinase subtype I  $\alpha$ , 8-Br-PET-cGMP is one of the best activators of isozyme I  $\beta$  and one of the best cyclic nucleotide muscle relaxants. BIOLOG also offers the corresponding PDE-stable structure Rp-8-Br-PET-cGMPS (Prod. No. [BLG-P007](#)) and Sp-8-Br-PET-cGMPS (Prod. No. [BLG-P008](#)).

*BLG-P003-50 (5 x 10  $\mu$ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.*

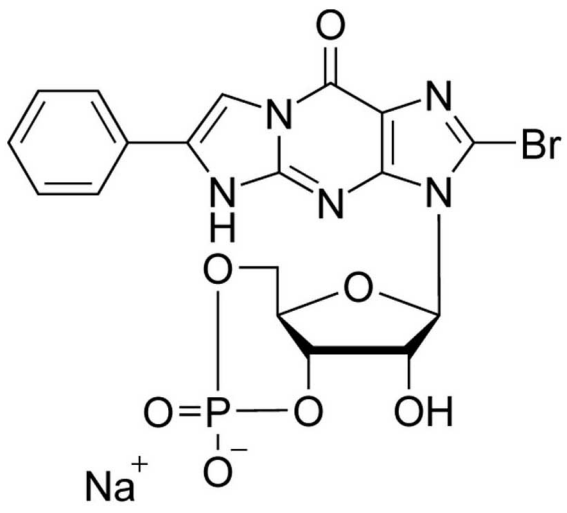
## Ordering Information

[Order Online »](#)

BLG-P003-10	10 $\mu$ mol
BLG-P003-50	5x10 $\mu$ 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**  $\beta$ -Phenyl-1,N2-etheno-8-bromoguanosine-3',5'-cyclic monophosphate . sodium salt

**CAS** 144510-04-3

**Couple Type** Modified nucleotides

**Formula**  $C_{18}H_{14}BrN_5O_7P \cdot Na$

**MW** 546.2

**Purity** > 98% HPLC

**Quantity** 10 $\mu$ mol $\approx$ 5.5mg

**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)