

Rp-8-Br-PET-cGMPS

Membrane-permeant and PDE-resistant PKG inhibitor

Competitive inhibitor of protein kinase G type I alpha and I beta and of cGMP-gated retinal type ion channels (cyclic GMP antagonist). Resistant against mammalian cyclic nucleotide-dependent phosphodiesterases, no metabolic side effects. More lipophilic and membrane permeant compared to Rp-8-pCPT-cGMPS (Prod. No. [BLG-C013](#)).

BLG-P007-05 (5 x 1 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

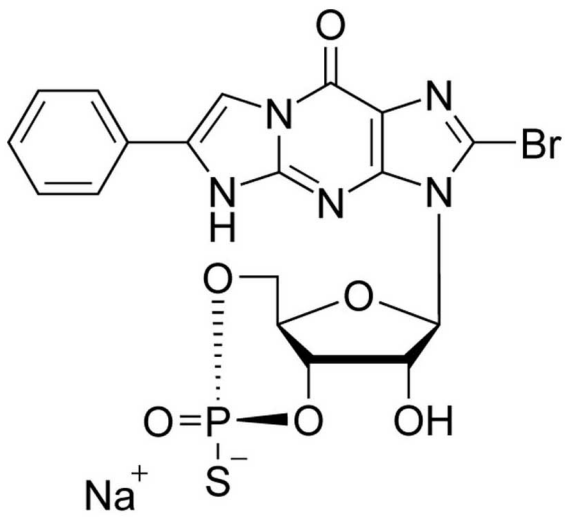
Ordering Information

[Order Online »](#)

BLG-P007-01	1µmol
BLG-P007-05	5x1µmol

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 β -Phenyl-1,N2-etheno-8-bromoguanosine-3',5'-cyclic monophosphorothioate . sodium salt, Rp-isomer

Appearance Lyophilized or crystallized sodium salt.

CAS 172806-20-1

Couple Type Modified nucleotides

Formula $C_{18}H_{14}BrN_5O_6PS \cdot Na$

MW 562.3

Purity > 99% HPLC

Quantity 1 μ mol \approx 0.6mg

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

Solubility Limited solubility 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

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com