

Sp-8-Br-PET-cGMPS

Membrane-permeant and PDE-resistant PKG agonist

Membrane-permeant inhibitor of the retinal type cGMP-gated ion channel (cGMP antagonist) but activator of cyclic GMP-dependent protein kinase I α and I β (cGMP agonist). Suitable for distinction between kinase and ion channel mediated effects. Resistant against mammalian cyclic nucleotide-dependent phosphodiesterase, no metabolic side effects. More lipophilic and permeant as compared to Sp-8-pCPT-cGMPS (Prod. No. [BLG-C014](#)).

BLG-P008-25 (5 x 5 μ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

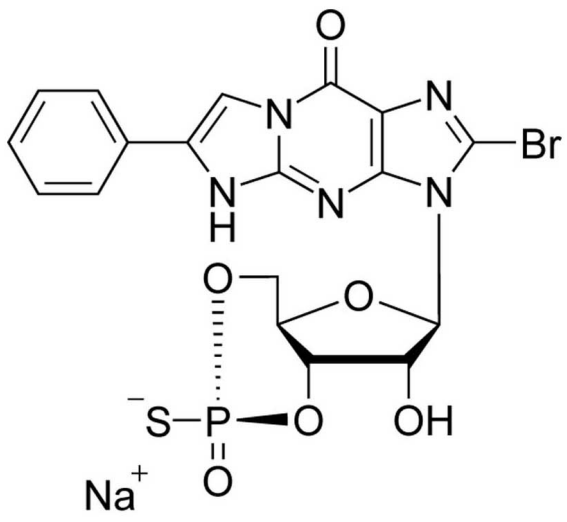
Ordering Information

[Order Online »](#)

BLG-P008-05	5 μ mol
BLG-P008-25	5x5 μ 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 β -Phenyl-1,N2-etheno-8-bromoguanosine-3',5'-cyclic monophosphorothioate . sodium salt, Sp-isomer

CAS 172806-21-2

Couple Type Activator, Modified nucleotides

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

MW 562.3

Purity > 99% HPLC

Quantity 5 μ mol \approx ~2.8mg

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