

Rp-8-pCTP-PET- cGMPS

Potent, PDE-resistant inhibitor of protein kinase G

Potent inhibitor of protein kinase G. Most probably retinal cGMP-gated ion channels inhibitor. Resistant against mammalian cyclic nucleotide-dependent phosphodiesterase.

BLG-C046-05 (5 x 1 μ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

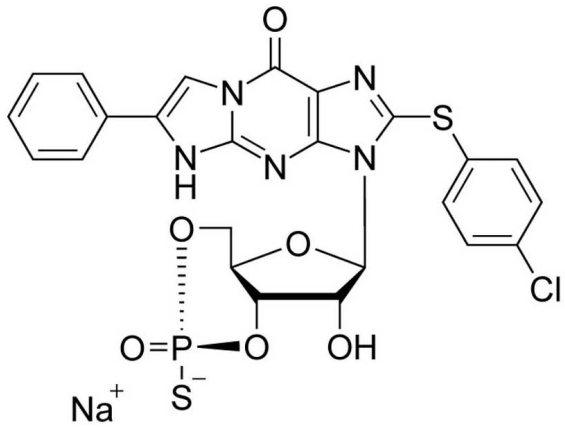
Ordering Information

[Order Online »](#)

BLG-C046-01	1 μ mol
BLG-C046-05	5x1 μ 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)-β-phenyl-1,N2-ethenoguanosine-3',5'-cyclic monophosphorothioate . sodium salt, Rp-isomer

CAS 1262749-62-1

Couple Type Modified nucleotides

Formula $C_{24}H_{18}ClN_5O_6PS_2 \cdot Na$

MW 626

Purity > 99% HPLC

Quantity 1μmol≈0.6mg

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