

PET-cGMP

Site-selective analogue of cyclic GMP

Potent membrane permeant activator of both isozymes I α and I β of cGMP-dependent protein kinase while only a poor stimulator of type II. Common agonists such as 8-Br-cGMP (Prod. No. [BLG-B004](#)) or 8-pCPT-cGMP (Prod. No. [BLG-C009](#)) mainly activate kinase subtype I α and type II.

BLG-P001-50 (5 x 10 μ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

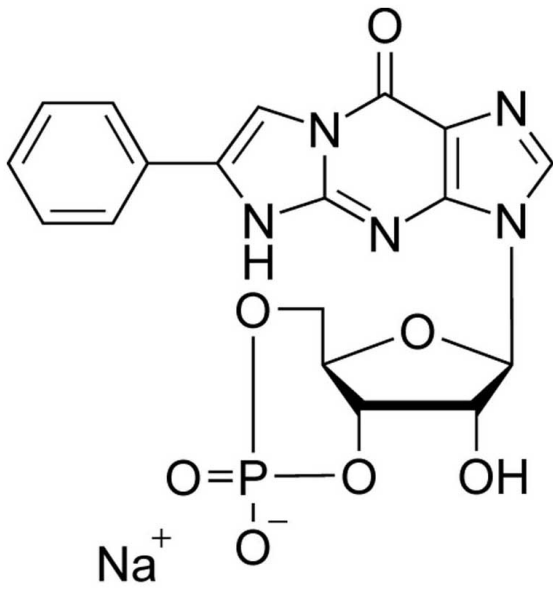
Ordering Information

[Order Online »](#)

BLG-P001-10	10 μ mol
BLG-P001-50	5x10 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-ethenoguanosine-3',5'-cyclic monophosphate . sodium salt

CAS 78080-27-0

Couple Type Activator, Modified nucleotides

Formula $C_{18}H_{15}N_5O_7P \cdot Na$

MW 467.3

Purity > 98% HPLC

Quantity 10 μ mol \approx 4.7mg

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