

# PET-cGMP-8-TMAmd-(EO)<sub>8</sub>-EAmdMT-8-cGMP-PET

## Polymer-linked dimeric (PLD) cGMP analogue

Polymer-linked dimeric (PLD) cGMP analogue in which two cGMP units are linked by a spacer with 8 ethylene oxide units via thioethers at the C8 positions. In addition,  $\beta$ -phenyl-1,N2-etheno (PET) groups have been attached to both guanine nucleobases.

Membrane-permeant cGMP dimer with enhanced PKG activation properties by addressing two binding sites simultaneously with a single molecule. Activation constants of PET-cGMP-8-TMAmd-(EO)<sub>8</sub>-EAmdMT-8-cGMP-PET in vitro on purified cGMP-dependent protein kinase (PKG) isozymes were measured to be in the range of < 1 nM (PKG1 $\alpha$  and PKG1 $\beta$ ) and ~35 nM (PKG2), respectively (BIOLOG in-house data).

*BLG-P054-05 (5 x 1  $\mu$ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.*

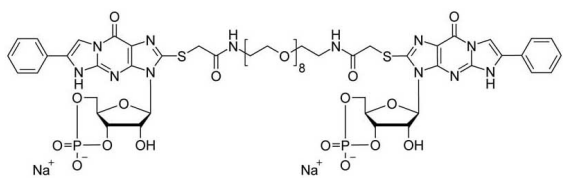
## Ordering Information

[Order Online »](#)

BLG-P054-01	1 $\mu$ mol
BLG-P054-05	5x1 $\mu$ 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-ethenoguanosine-3',5'-cyclic monophosphate-[8-thiomethylamido-(octaethoxy)-ethylamidomethylthio-8]- $\beta$ -phenyl-1,N2- ethenoguanosine-3',5'-cyclic monophosphate . sodium salt

**Couple Type** Modified nucleotides

**Formula**  $C_{58}H_{70}N_{12}O_{24}P_2S_2 \cdot Na_2$

**MW** 1491.3

**Purity** > 98% HPLC

**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)