

PET-cGMP-8-TMAmd- (PEG pd 2000)- AmdMT-8-cGMP-PET

Polymer-linked dimeric (PLD) cGMP analogue

Polymer-linked dimeric (PLD) cGMP analogue in which two cGMP units are linked by a polydisperse 2000 Dalton spacer on basis of polyethylene glycol (PEG) via thio-functions at the C8 positions. In addition, β -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-(PEG pd 2000)-AmdMT-8-cGMP-PET in vitro on purified cGMP-dependent protein kinase (PKG) isozymes were measured to be in the range of < 1 nM (PKG1 α and PKG1 β) and ~80 nM (PKG2), respectively (BIOLOG in-house data). PET-cGMP-8-TMAmd-(PEG pd 2000)-AmdMT-8-cGMP-PET decreased cell viability in the PKG2 expressing colon cancer cell line Caco-2, but not in the PKG2 expressing colon cancer cell lines HCT 116 and HT-29.

*BLG-P059-05 (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-P059-01	1 μ mol
BLG-P059-05	5x1 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-[8-thiomethylamido-(PEG pd 2000)-amidomethylthio-8]- β -phenyl-1,N2-ethenoguanosine-3',5'-cyclic monophosphate . sodium salt

Couple Type Modified nucleotides

Formula $\sim C_{130}H_{216}N_{12}O_{60}P_2S_2$ (free acid)

MW ~ 3033

Purity > 95% HPLC (est.)

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