

# cGMP-8-T-(EO)<sub>5</sub>-ET-8-cGMP

## Polymer-linked dimeric (PLD) cGMP analogue

Polymer-linked dimeric (PLD) cGMP analogue in which two cGMP units are linked by a spacer with 5 ethylene oxide units via thioethers at the C8 positions. Membrane-permeant cGMP dimer with enhanced PKG activation properties by addressing two binding sites simultaneously with a single molecule.

Activation constants of cGMP-8-T-(EO)<sub>5</sub>-ET-8-cGMP in vitro on purified cGMP-dependent protein kinase (PKG) isozymes were measured to be in the range of < 1 nM (PKG1 $\beta$  and PKG2) and < 0.8 nM (PKG1 $\alpha$ ), respectively. The compound significantly reduced proliferation and migration combined with increased cell death in two melanoma cell lines. In the colon cancer cell line HT-29, treatment with cGMP-8-T-(EO)<sub>5</sub>-ET-8-cGMP led to a significant reduction of cell viability.

*BLG-G031-05 (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-G031-01	1 $\mu$ mol
BLG-G031-05	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** Guanosine-3',5'-cyclic monophosphate-[8-thio-(pentaethoxy)-ethylthio-8]-guanosine-3',5'-cyclic monophosphate . sodium salt

**CAS** 2156612-58-5

**Couple Type** Modified nucleotide

**Formula**  $C_{32}H_{44}N_{10}O_{19}P_2S_2 \cdot Na_2$

**MW** 1044.8

**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)