

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

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 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-TMAmd-(PEG pd 2000)-AmdMT-8-cGMP in vitro on purified cGMP-dependent protein kinase (PKG) isozymes were measured to be in the range of 7.7 ± 1.3 nM (PKG1 α), 21 ± 2 nM (PKG1 β), and 2.1 ± 0.6 nM (PKG2), respectively. The compound significantly reduced proliferation, migration and increased cell death in two melanoma cell lines. In the colon cancer cell lines Caco-2 and HCT 116, treatment with cGMP-8-TMAmd-(PEG pd 2000)-AmdMT-8-cGMP led to a significant reduction of cell viability.

BLG-G043-05 (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-G043-01	1 μ mol
BLG-G043-05	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 Guanosine-3',5'-cyclic monophosphate-[8-thiomethylamido-(PEG pd 2000)-amidomethylthio-8]-guanosine-3',5'-cyclic monophosphate . sodium salt

Couple Type Modified nucleotide

Formula ~ C₁₁₄H₂₀₈N₁₂O₆₀P₂S₂ (free acid)

MW ~ 2833 (free acid)

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