

# Di-MANT-c-diGMP

## Fluorescent analogue of c-diGMP

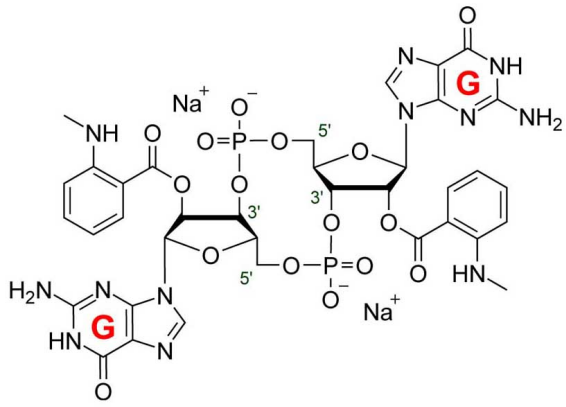
Di-MANT-c-diGMP is a fluorescent analogue of the natural signalling molecule c-diGMP (Prod. No. BLG-C057). The MANT fluorophore ( $\lambda_{\text{exc}} = 355 \text{ nm}$ ,  $\lambda_{\text{em}} = 448 \text{ nm}$ ) has a certain sensitivity for its environment and can change its spectral properties upon binding.

*BLG-D101-005 (5 x 0.1  $\mu\text{mol}$  pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.*

### Ordering Information [Order Online »](#)

BLG-D101-001	0.1 $\mu\text{mol}$
BLG-D101-005	5x0.1 $\mu\text{mol}$

### Manuals, SDS & CofA [View Online »](#)



## Handling & Storage

Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

Alternative Name	2',2"-O-(Di-N'-methylantraniloyl)-cyclic diguanosine monophosphate . sodium salt
Couple Type	Modified nucleotides
Formula	$C_{36}H_{38}N_{12}O_{16}P_2$ (free acid)
MW	956.7 (free acid)
Purity	> 95% HPLC
Quantity	0.1µmol≈~0.1mg
Solubility	Very soluble in water.
Technical Info / Product Notes	For the Original Manufacturer's data sheet please <a href="#">click here</a> .



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)