

# MANT-GppNHp

## Fluorescent form of the hydrolysis-resistant GppNHp

Fluorescent analogue with  $\lambda_{\text{exc}}$  355 nm and  $\lambda_{\text{em}}$  448 nm, useful for research into GTP-dependent receptor proteins. The MANT fluorophore has a certain sensitivity for its environment and can change its spectral properties upon binding. Potent inhibitor of adenylate cyclase.

*BLG-M047-05 (5 x 1  $\mu\text{mol}$  pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.*

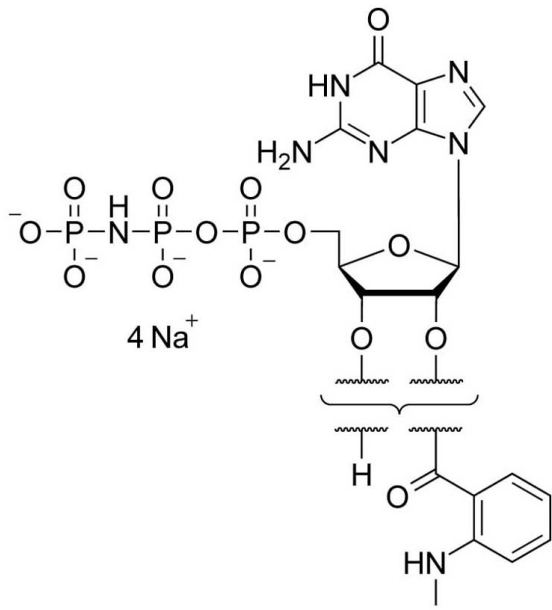
### Ordering Information

[Order Online »](#)

BLG-M047-01	1 $\mu\text{mol}$
BLG-M047-05	5x1 $\mu\text{mol}$

### Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

<b>Use/Stability</b>	Compound slowly decomposes at ambient temperature.
<b>Handling</b>	Protect from light.
<b>Long Term Storage</b>	-80°C
<b>Shipping</b>	Dry Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	2'/3'-O-(N'-Methylantraniloyl)guanosine-5'-O-[(β,γ)-imido]triphosphate . tetrasodium salt
<b>Appearance</b>	10 mM aqueous solution
<b>CAS</b>	148821-01-6
<b>Couple Type</b>	Modified nucleotides
<b>Formula</b>	$C_{18}H_{24}N_7O_{14}P_3$ (free acid)
<b>MW</b>	655.4
<b>Purity</b>	> 90% HPLC
<b>Quantity</b>	1 μmol ≈ 0.7mg
<b>Technical Info / Product Notes</b>	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)