

# MANT-GMP

## Fluorescent analogue of 5'-GMP

Fluorescent metabolite of MANT-cGMP and probe for 5'-GMP-binding receptors. Analog of guanosine-5'-monophosphate, useful for research into 5'-GMP-dependent receptor proteins.

*BLG-M042-50 (5 x 10  $\mu$ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.*

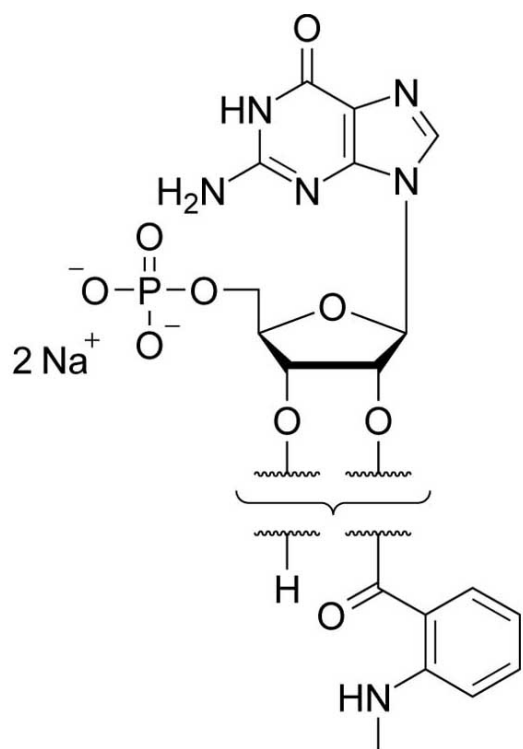
### Ordering Information

[Order Online »](#)

<b>BLG-M042-10</b>	10 $\mu$ mol
<b>BLG-M042-50</b>	5x10 $\mu$ 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'-/3'-O-(N'-Methylantraniloyl)guanosine-5'-O-monophosphate . disodium salt
CAS	85287-54-3
Couple Type	Modified nucleotides
Formula	$C_{18}H_{21}N_6O_9P$ (free acid)
MW	496.4 (free acid)
Purity	> 97% HPLC for mixture of isomers
Quantity	10µmol≈5mg
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-fr@enzolifesciences.com](mailto:info-fr@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)