

MANT-c-diGMP

Fluorescent analogue of c-diGMP

Fluorescent analogue of the bacterial second messenger c-diGMP (Cat No. BLG-C057). The MANT fluorophore (λ_{exc} 355 nm; λ_{em} 448 nm) has a certain sensitivity for its environment and can change its spectral properties upon binding. Compare: Hiratsuka, *Biochim. Biophys. Acta*, **742**, 496 - 503 (1983).

BLG-M102-005 (5 x 0.1 μ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

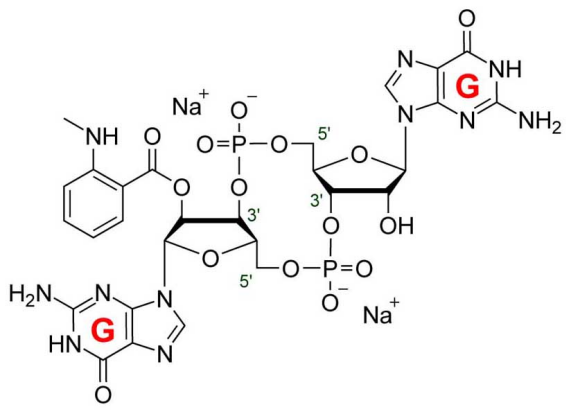
Ordering Information

[Order Online »](#)

BLG-M102-001	0.1 μ mol
BLG-M102-005	5x0.1 μ mol

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	Light-sensitive.
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	2'-O-(N'-Methylantraniloyl)-cyclic diguanosine monophosphate . sodium salt
CAS	1384529-33-2
Formula	$C_{28}H_{31}N_{11}O_{15}P_2$ (free acid)
MW	823.6 (free acid)
Purity	> 95% HPLC
Quantity	0.1 μmol ≈ 82 μg
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