

MANT-c-diAMP

Fluorescent analogue of c-diAMP

*BLG-M103-005 (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-M103-001	0.1 μ mol
BLG-M103-005	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 diadenosine monophosphate . sodium salt
Formula	$C_{28}H_{31}N_{11}O_{13}P_2$ (free acid)
MW	791.6 (free acid)
Purity	> 95% HPLC
Quantity	0.1 μmol ≈ 79 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](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)