

MANT-cCMP

Fluorescent, membrane-permeant cyclic
CMP analogue for PDE studies

Membrane-permeant, fluorescent analogue with $\lambda_{\text{exc}} = 350 \text{ nm}$ and $\lambda_{\text{em}} = 441 \text{ nm}$, potentially suitable for phosphodiesterase studies. The compound is expected to show a decrease in fluorescence when cleaved by PDE.

*BLG-M053-25 (5 x 5 μmol pack size) is not sold in the U.S. or Canada.
Please [contact us](#) for available options.*

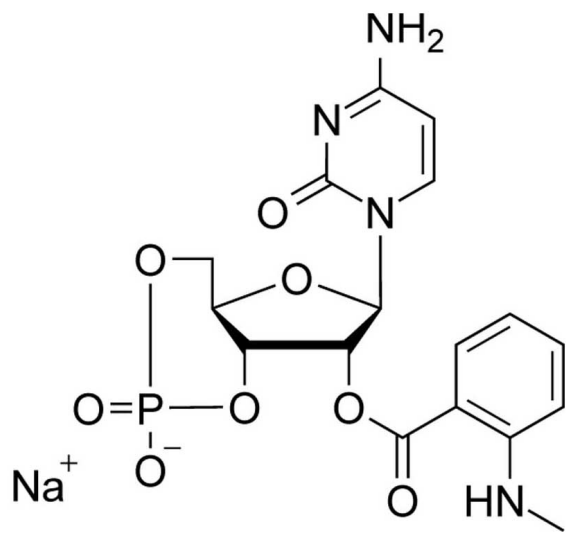
Ordering Information

[Order Online »](#)

BLG-M053-05	5 μmol
BLG-M053-25	5x5 μmol

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'-O-(N'-Methylantraniloyl)cytidine-3',5'-cyclic monophosphate . sodium salt
CAS	1477570-24-3
Couple Type	Modified nucleotides
Formula	$C_{17}H_{18}N_4O_8P \cdot Na$
MW	460.3
Purity	> 98% HPLC
Quantity	5µmol≈2.3mg
Solubility	Soluble in water.

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