

Mant-cAMP

Fluorescent, PKA-inactive analogue of cyclic AMP

Fluorescent, membrane-permeant cyclic AMP analogue, suitable for phosphodiesterase studies.

The compound shows a decrease in fluorescence when cleaved by PDE.

BLG-M008-50 (5 x 10 μ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

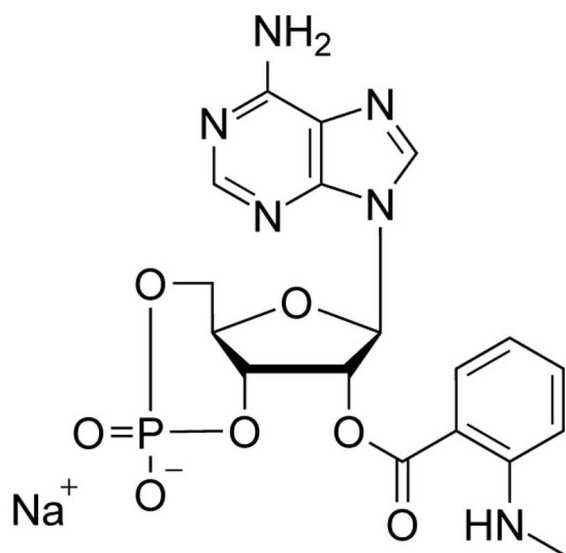
Ordering Information

[Order Online »](#)

BLG-M008-10	10 μ mol
BLG-M008-50	5x10 μ 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'-O-(N'-Methylantraniloyl)adenosine-3',5'-cyclic monophosphate . sodium salt
CAS	83707-14-6
Couple Type	Modified nucleotides
Formula	$C_{18}H_{18}N_6O_7P \cdot Na$
MW	484.4
Purity	> 98% HPLC
Quantity	10µmol≈4.8mg
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-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-de@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-uk@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com