

6-MAH-cAMP

Fluorescent cyclic AMP analogue

Fluorescent cyclic AMP analogue. Excitation 333nm, Emission 437nm.

*BLG-M017-25 (5 x 5 μ mol pack size) is not sold in the U.S. or Canada.
Please [contact us](#) for available options.*

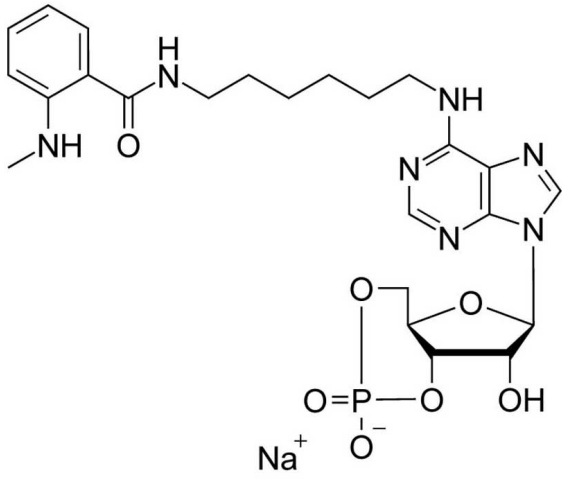
Ordering Information

[Order Online »](#)

BLG-M017-05	5 μ mol
BLG-M017-25	5x5 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	N6-(6-[N ¹ -Methylanthraniloyl]aminohexyl)adenosine-3',5'-cyclic monophosphate . sodium salt
CAS	723313-27-7
Couple Type	Modified nucleotides
Formula	C ₂₄ H ₃₁ N ₇ O ₇ P . Na
MW	583.5
Purity	> 98% HPLC
Quantity	5µmol≈2.9mg
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