

MANT-AMP

Fluorescent analogue of 5'-AMP

Fluorescent metabolite of MANT-cAMP and probe for 5'-AMP-binding receptors. Analog of adenosine-5'-monophosphate, useful for research into 5'-AMP-dependent receptor proteins.

BLG-M037-50 (5 x 10 μ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

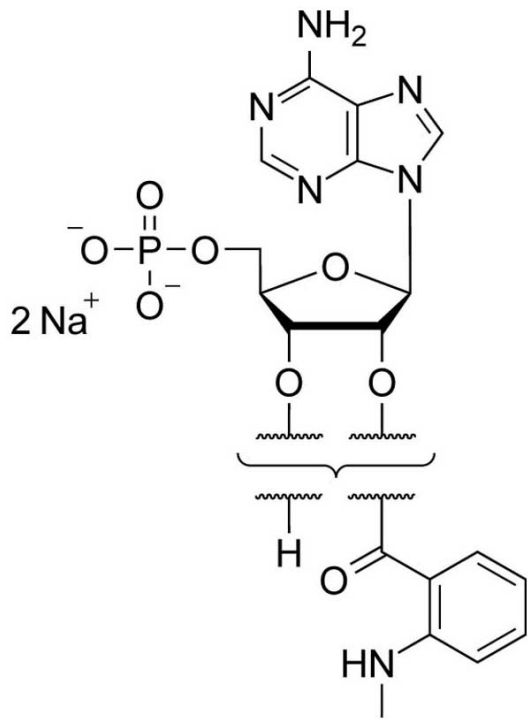
Ordering Information

[Order Online »](#)

BLG-M037-10	10 μ mol
BLG-M037-50	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'-/3'-O-(N'-Methylantraniloyl)adenosine-5'-O-monophosphate . disodium salt

CAS 85287-53-2

Couple Type Modified nucleotides

Formula $C_{18}H_{21}N_6O_8P$ (free acid)

MW 480.4 (free acid)

Purity > 97% HPLC for mixture of isomers

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-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