

MANT- Ap_4 -MANT-A

BLG-M131-05 (5 x 1 μmol pack size) is not sold in the U.S. or Canada.
Please [contact us](#) for available options.

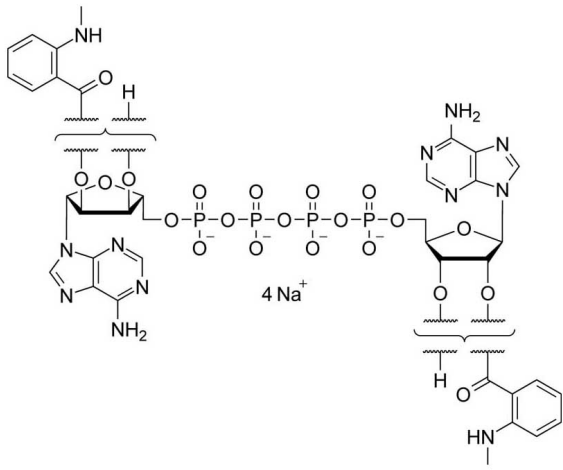
Ordering Information

[Order Online »](#)

BLG-M131-01	1 μmol
BLG-M131-05	5x1 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	P1-(5'-[2'-/3'-O-(N'-Methylanthraniloyl)adenosyl])-P4-(5'-[2'-/3'-O-(N'-methylanthraniloyl)adenosyl])-tetraphosphate . tetrasodium salt, MANT-Apppp-MANT-A
Couple Type	Modified nucleotides
Formula	$C_{36}H_{42}N_{12}O_{21}P_4$ (free acid)
MW	1102.7 (free acid)
Purity	> 95% HPLC
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