

MANT-5'-IMP

Fluorescent analogue of 5'-IMP

Fluorescent analogue of inosine-5'-monophosphate (λ_{exc} 355 nm, λ_{em} 448 nm). The MANT fluorophore has a certain sensitivity for its environment and can change its spectral properties upon binding.

BLG-M072-50 (5 x 10 μ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

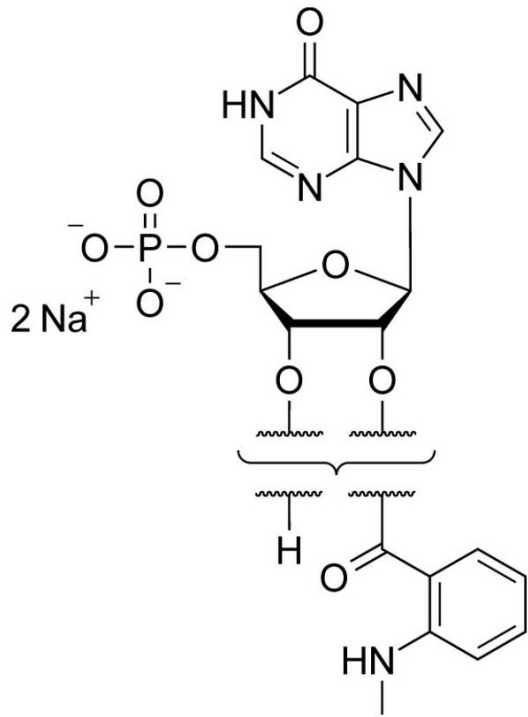
Ordering Information

[Order Online »](#)

BLG-M072-10	10 μ mol
BLG-M072-50	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)inosine-5'-O-monophosphate . disodium salt
Couple Type	Modified nucleotides
Formula	$C_{18}H_{20}N_5O_9P$ (free acid)
MW	481.4
Purity	> 97% HPLC for mixture of isomers



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