

8-Br-2'-O-Me-cGMP

Epac-negative control

*BLG-B037-25 (5 x 5 μ mol pack size) is not sold in the U.S. or Canada.
Please [contact us](#) for available options.*

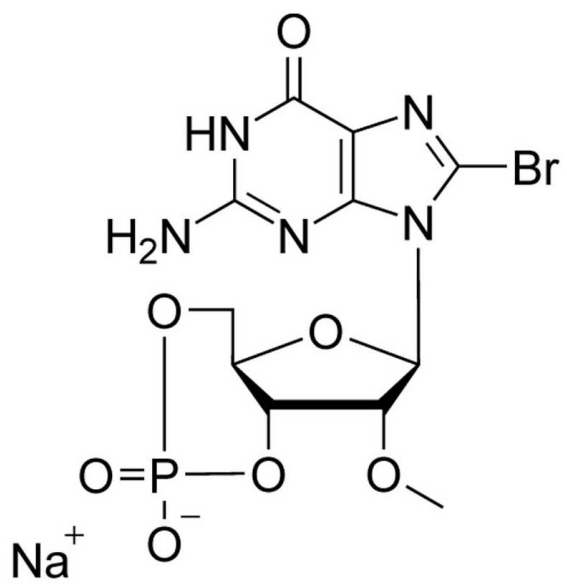
Ordering Information

[Order Online »](#)

BLG-B037-05	5 μ mol
BLG-B037-25	5x5 μ mol

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 8-Bromo-2'-O-methylguanosine-3',5'-cyclic monophosphate . sodium salt

Couple Type Modified nucleotides

Formula $C_{11}H_{12}BrN_5O_7P \cdot Na$

MW 460.1

Purity > 98% HPLC

Quantity 5µmol≈1.9mg

Technical Info / Product Notes **Legal information:** the reagent is subject to patent application. It is sold under limited, non transferable licence for research purposes only, to the exclusion of any commercial use.



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