

8-Br-c-diGMP

Mono-functionalized analogue of c-diGMP

Mono-functionalized analogue of the bacterial second messenger c-diGMP (Cat. No. BLG-C057), useful for binding studies and precursor for further modifications with spacers or labels.

BLG-B099-005 (5 x 0.1 μ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

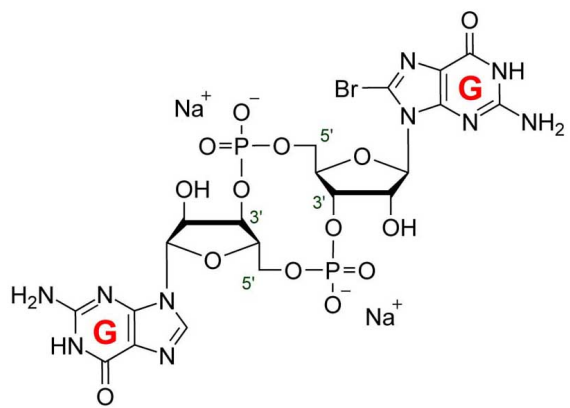
Ordering Information

[Order Online »](#)

BLG-B099-001	0.1 μ mol
BLG-B099-005	5x0.1 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-cyclic diguanosine monophosphate (8-Br-c-diGMP) . sodium salt

Couple Type Modified nucleotides

Formula $C_{20}H_{23}BrN_{10}O_{14}P_2$ (free acid)

MW 769.3 (free acid)

Purity > 95% HPLC

Quantity 0.1 μmol ≈ 77 μg

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