

2'-Biotin-16-c-diGMP

Biotinylated analogue of the bacterial second messenger c-diGMP

Analogue of c-diGMP with biotin conjugate most probably suitable as tracer in immunoassays.

BLG-B098-005 (5 x 0.1 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

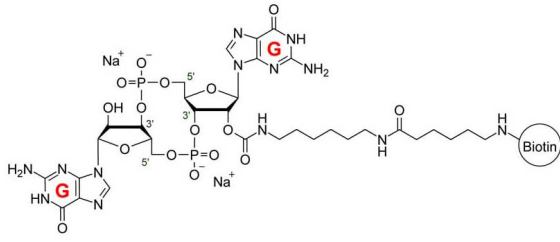
Ordering Information

[Order Online »](#)

BLG-B098-001	0.1µmol
BLG-B098-005	5x0.1µmol

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 2'-O-(6-(6-[Biotinyl]aminohexanoyl)aminohexylcarbamoyl)-cyclic diguanosine monophosphate . sodium salt

Couple Type Modified nucleotides

Formula $C_{43}H_{63}N_{15}O_{18}P_2S$ (free acid)

MW 1172.1 (free acid)

Purity > 95% HPLC

Quantity 0.1 µmol ≈ 0.1 mg

Solubility Soluble in water or in aqueous buffers (>3.5mM).

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