

c-(Ap-8-Biotin-11-Gp)

Biotinylated analogue of the bacterial second messenger cGAMP

c-(Ap-8-[Biotin]-AET-Gp) is a biotinylated analogue of the cyclic dinucleotide c-(ApGp) (*aka* cGAMP, Prod. No. BLG-C117) in which the biotin label is attached to position 8 of the guanine nucleobase via a 11 atom spacer.

BLG-C157-005 (5 x 0.1 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

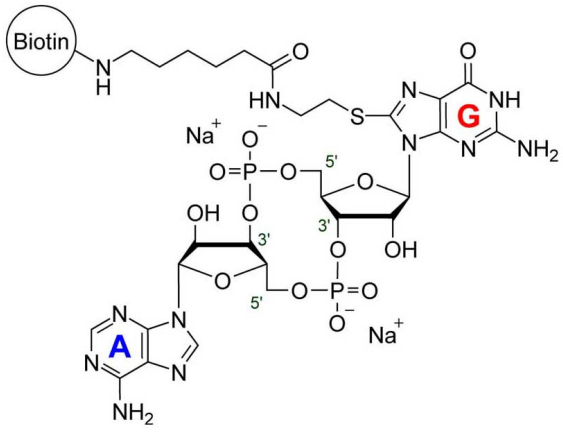
Ordering Information

[Order Online »](#)

BLG-C157-001	0.1µmol
BLG-C157-005	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 Cyclic (adenosine monophosphate-8-(2-(6-[biotinyl]aminohexanoyl)aminoethylthio)guanosine monophosphate) . sodium salt

Formula $C_{38}H_{54}N_{14}O_{16}P_2S_2$ (free acid)

MW 1089.0 (free acid)

Purity > 95% HPLC

Solubility Soluble in water and aqueous buffers (>1.5mM).

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