

c[8-Biotin-11- G(2',5')pA(3',5')p]

Biotinylated analogue of the metazoan
second messenger 2',3'-cGAMP

BLG-C176-005 (5 x 0.1 µmol pack size) is not sold in the U.S. or Canada.

Please [contact us](#) for available options.

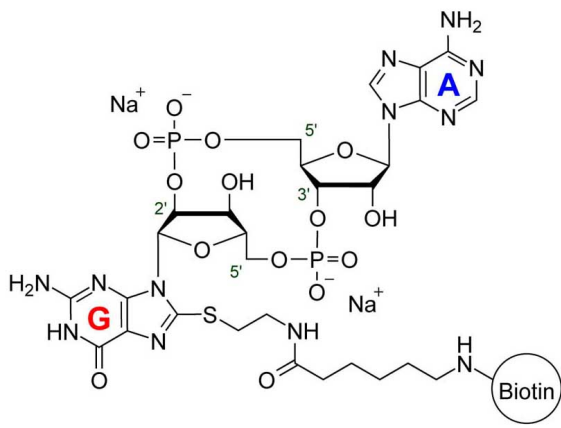
Ordering Information

[Order Online »](#)

BLG-C176-001	0.1µmol
BLG-C176-005	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 (8-(2-(6-[biotinyl]aminohexanoyl)aminoethylthio)guanosine-(2'→5')-monophosphate-adenosine-(3→5')-monophosphate) . sodium salt

Formula $C_{38}H_{54}N_{14}O_{16}P_2S_2$

MW 1089.0 (free acid)

Purity > 95% HPLC

Quantity ~0.1 mg

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