

# c[G(2',5')p-2'-[Biotin]- AHC-A(3',5')p]

BLG-C197-005 (5 x 0.1  $\mu$ mol pack size) is not sold in the U.S. or Canada.

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

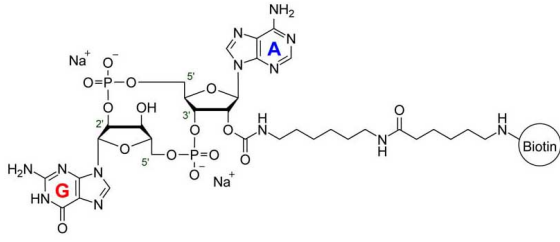
## Ordering Information

[Order Online »](#)

BLG-C197-001	0.1 $\mu$ mol
BLG-C197-005	5x0.1 $\mu$ 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** Cyclic(guanosine-(2'→5')-monophosphate-2'-O-(6-[biotinyl]aminohexylcarbamoyl)adenosine-(3'→5')-monophosphate) . sodium salt, c[G(2',5')p-2'-[Biotin]-AHC-A(3',5')p]

**Formula**  $C_{43}H_{63}N_{15}O_{17}P_2S$  (free acid)

**MW** 1156.1

**Purity** > 95% HPLC

**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](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)