

c[3'-Biotin-16- G(2',5')pA(3',5')p]

BLG-C196-005 (5 x 0.1 μ mol pack size) is not sold in the U.S. or Canada.
Please [contact us](#) for available options.

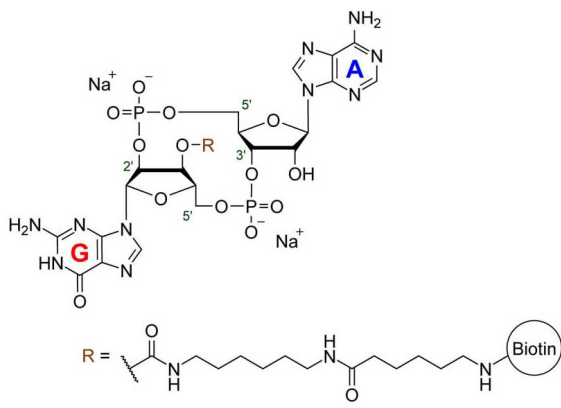
Ordering Information

[Order Online »](#)

BLG-C196-001	0.1 μ mol
BLG-C196-005	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 Cyclic (3'-O-(6-(6-[biotinyl]amino)hexanoyl)amino)hexylcarbamoyl)guanosine-(2'→5')-monophosphate-adenosine-(3'→5')-monophosphate) . sodium salt, c[3'-[Biotin]-AHC-G(2',5')pA(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

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