

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

Metazoan cyclic dinucleotide second messenger

c[G(2',5')pA(3',5')p] is a non-canonical cyclic dinucleotide in which a 5'-GMP unit is connected with a 5'-AMP unit via a 2'-5' and a 3'-5' phosphodiester bond to form a cyclic structure.

BLG-C161-05 (5 x 1 μ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

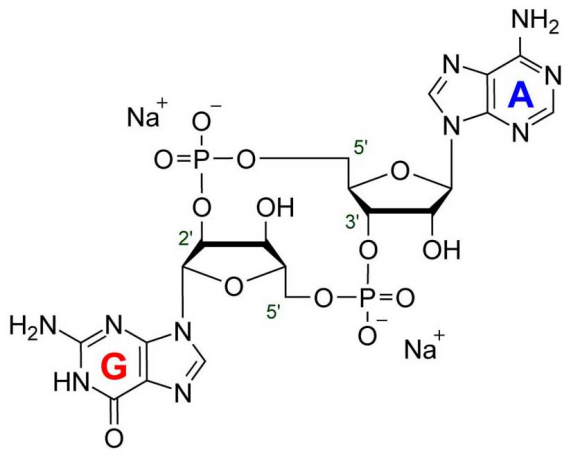
Ordering Information

[Order Online »](#)

BLG-C161-01	1 μ mol
BLG-C161-05	5x1 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 (guanosine-(2'→5')-monophosphate-adenosine-(3'→5')-monophosphate) . sodium salt

CAS 1441190-66-4

Formula $C_{20}H_{24}N_{10}O_{13}P_2$ (free acid)

MW 674.4 (free acid)

Purity > 98% HPLC

Quantity 0.5µmol≈335µg

Solubility Soluble in water and in aqueous buffers (>>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