

# c[A(2',5')pG(2',5')p]

BLG-C210-05 (5 x 1  $\mu$ mol pack size) is not sold in the U.S. or Canada.  
Please [contact us](#) for available options.

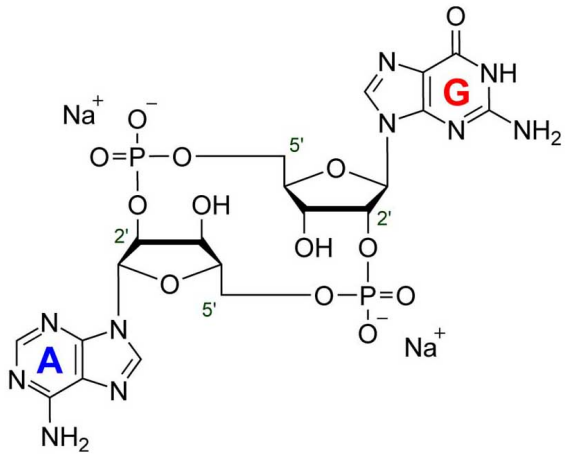
## Ordering Information

[Order Online »](#)

BLG-C210-01	1 $\mu$ mol
BLG-C210-05	5x1 $\mu$ mol

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

<b>Use/Stability</b>	c[A(2',5')pG(2'5')p] has sufficient stability at room temperature, and does not need special care during handling or shipment. It is recommended that the compound should be stored in the freezer, for longer storage periods.
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	Cyclic (adenosine-(2'→5')-monophosphate-guanosine-(2'→5')-monophosphate) . sodium salt, 2'2'-cGAMP, 2',5'-2',5'-cGAMP
<b>CAS</b>	1465774-27-9
<b>Formula</b>	$C_{20}H_{24}N_{10}O_{13}P_2$ (free acid)
<b>MW</b>	674.4 (free acid)
<b>Purity</b>	> 98% HPLC
<b>Technical Info / Product Notes</b>	For the Original Manufacturer's data sheet please <a href="#">click here</a> .



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)