

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

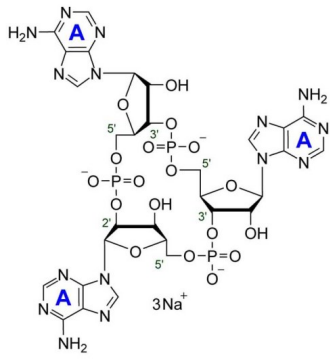
c[A(2',5')pA(3',5')pA(3',5')p] was identified as a bacterial signalling nucleotide involved in CBASS (cyclic oligonucleotide-based antiphage signalling system) immunity which protects prokaryotes from phage infection.

*BLG-C372-025 (5 x 0.5 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.*

## Ordering Information [Order Online »](#)

BLG-C372-005	0.5µmol
BLG-C372-025	5x0.5µmol

## Manuals, SDS & CofA [View Online »](#)



## Handling & Storage

Long Term Storage -20°C

Shipping Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	Cyclic (adenosine-(2'→5')-monophosphate-adenosine-(3'→5')-monophosphate-adenosine-(3'→5')-monophosphate) . sodium salt, 2'3'3'-cAAA, 2'-5',3'-5',3'-5'-c-AMP-AMP-AMP
<b>Appearance</b>	Solid
<b>CAS</b>	153736-11-9
<b>Couple Type</b>	Modified nucleotides
<b>Formula</b>	$C_{30}H_{36}N_{15}O_{18}P_3$ (free acid)
<b>MW</b>	987.6
<b>Purity</b>	> 95% HPLC
<b>Quantity</b>	0.5µmol are equivalent to ~0.5 mg



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)