

# c[A(3',5')pA(3',5')pG(3',5')p] / cAAG

## Bacterial cyclic trinucleotide

c[A(3',5')pA(3',5')pG(3',5')p] is a cyclic trinucleotide that was found to be biosynthesized by CD-NTase038 from the *Enterobacter cloacae* complex (strain UCI 50) (WP\_032676400, also named EcCdnD02). While c[A(3',5')pA(3',5')pG(3',5')p] is not recognized by STING, it binds to the mammalian cyclic dinucleotide sensor RECON (REductase COntrolling NF-κB) (all data according to Whiteley et al., *Nature*, 567, 194 - 199 (2019)).

*BLG-C361-05 (5 x 1 μmol pack size) is not sold in the U.S. or Canada.  
Please [contact us](#) for available options.*

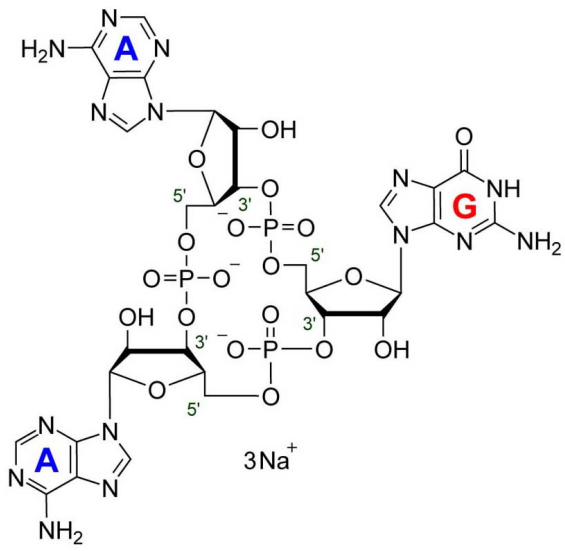
## Ordering Information

[Order Online »](#)

BLG-C361-01	1 μmol
BLG-C361-05	5x1 μmol

## 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 (adenosine-(3'→5')-monophosphate-adenosine-(3'→5')-monophosphate-guanosine-(3'→5')-monophosphate) . sodium salt, Cyclic AMP-AMP-GMP, 3'3'3'-cAAG

**CAS** 2365165-54-2

**Couple Type** Modified nucleotides

**Formula**  $C_{30}H_{36}N_{15}O_{19}P_3$  (free acid)

**MW** 1003.6 (free acid)

**Purity** > 95% HPLC

**Quantity** 0.5µmol is equivalent to ~0.5 mg

**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)