

# c-triAMP

## Cyclic oligoadenylate involved in Type III CRISPR-Cas systems in prokaryotes

c-triAMP is a cyclic oligoadenylate involved in Type III CRISPR-Cas systems in prokaryotes. Upon recognition and binding of invasive target RNA, the Cas10 subunit of the type III interference complex generates cyclic oligoadenylates which in turn act as allosteric activators of Csm6 ribonucleases that degrade invader-derived RNA transcripts. c-triAMP was found to be the predominant reaction product in *Streptococcus thermophilus* while only c-hexaAMP activates the Csm6 ribonuclease of this organism (Kazlauskienė et al. (2017)).

BLG-C362-05 (5 x 1 µmol pack size) is not sold in the U.S. or Canada.  
Please [contact us](#) for available options.

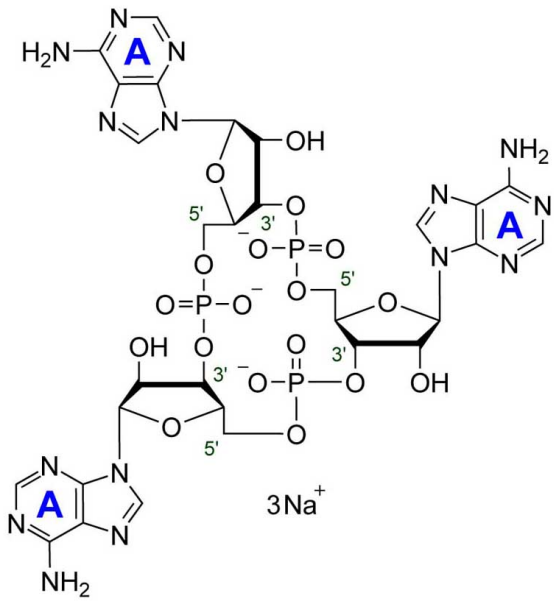
### Ordering Information

[Order Online »](#)

BLG-C362-01	1µmol
BLG-C362-05	5x1µmol

### 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 triadenosine monophosphate . sodium salt, Cyclic triadenylate, c-A3, cOA3

**Formula**  $C_{30}H_{36}N_{15}O_{18}P_3$  (free acid)

**MW** 987.6 (free acid)

**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](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)