

# c-hexaAMP

c-hexaAMP is a cyclic nucleotide in which six 5'AMP units are interconnected via 3'-5' phosphodiester bonds to form a cyclic structure.

*BLG-C332-025 (5 x 0.5  $\mu$ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.*

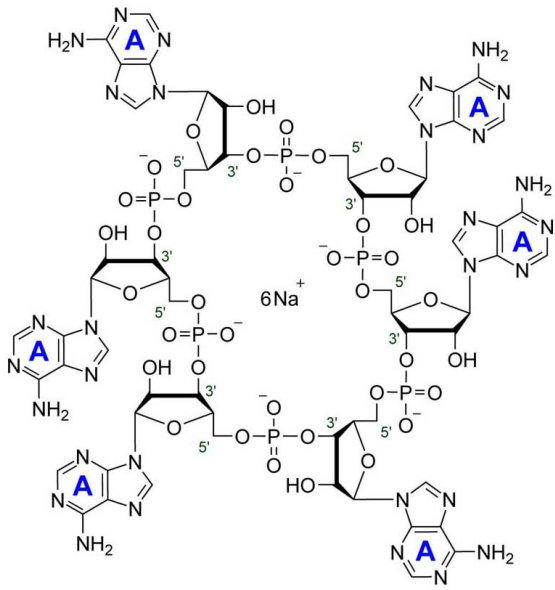
## Ordering Information

[Order Online »](#)

|              |                 |
|--------------|-----------------|
| BLG-C332-005 | 0.5 $\mu$ mol   |
| BLG-C332-025 | 5x0.5 $\mu$ mol |

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

|                          |  |
|--------------------------|--|
| <b>Use/Stability</b>     | c-hexaAMP has sufficient stability at room temperature and does not need special care during handling or shipment. Nevertheless, we recommend that the compound should be stored in the freezer, for longer storage periods preferably in freeze-dried form. |
| <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 hexaadenosine monophosphate . sodium salt, Cyclic hexaadenylate, c-A6  |
| <b>Appearance</b>         | Crystallized or lyophilized sodium salt   |
| <b>CAS</b>                | 232933-63-0   |
| <b>Endotoxin Content</b>  | Has not been tested for endotoxins.   |
| <b>Excitation Maximum</b> | 259 nm  |
| <b>Formula</b>            | $C_{60}H_{72}N_{30}O_{36}P_6$ (free acid)   |
| <b>MW</b>                 | 1975.3 (free acid)  |
| <b>Purity</b>             | > 95% HPLC  |
| <b>Solubility</b>         | Soluble in water and aqueous buffers. Please rinse tube walls carefully and preferably use ultrasonic or vortex to achieve total and uniform mixing. When opening the tube please make sure that no substance is lost within the cap. |

**Technical Info / Product Notes** Cyclic oligoadenylates such as c-hexaAMP were found to be novel bacterial second messengers involved in the Type 3 CRISPR-CAS associated detection and degradation of invasive genetic elements in many prokaryotes. Upon recognition and binding of invasive target RNA, the Cas10 subunit of the type 3 interference complex generates cyclic oligoadenylates which in turn act as allosteric activators of Csm6 ribonucleases that degrade invader-derived RNA transcripts. It is suggested that the size of cyclic oligoadenylate depends on the type 3 CRISPR-CAS system present, with c-hexaAMP being the predominant signaling molecule in *Enterococcus italicus* and *Streptococcus thermophilus* (all data according to Niewoehner et al., (2017) and Kazlauskienė et al., (2017)).



ENZO LIFE SCIENCES  
INC.  
Phone: 800.942.0430  
info-  
usa@enzolifesciences.com

European Sales Office  
ENZO LIFE SCIENCES  
(E.L.S.) AG  
Phone: +41 71 930 8389  
info-  
eu@enzolifesciences.com

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
info-  
be@enzolifesciences.com

France  
Phone: +33 472 440 633  
info-  
fr@enzolifesciences.com

Germany  
Phone: +49 621 6900 328  
info-  
de@enzolifesciences.com

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 823900  
info-  
uk@enzolifesciences.com