

# 2',2'',2''',2''''-Tetra-c-tetradAMP

Analogue of c-tetraAMP containing deoxy modifications in all ribose 2'-positions

Cyclic oligoadenylates such as c-tetraAMP (Biolog Cat. No. C 335) were found to be novel bacterial second messengers involved in the Type III CRISPR-Cas-associated detection and degradation of invasive genetic elements in many prokaryotes. 2',2'',2''',2''''-Tetra-c-tetradAMP is an analogue of c-tetraAMP containing deoxy modifications in all ribose 2'-positions, which are expected to make the compound resistant to degradation by nucleases.

*BLG-T070-025 (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-T070-005	0.5µmol
BLG-T070-025	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

**Alternative Name** 2'-, 2'', 2''', 2''''-Tetradeoxy-cyclic tetraadenosine monophosphate . sodium salt

**CAS** 125682-17-9

**Couple Type** Modified nucleotides

**Formula**  $C_{40}H_{48}N_{20}O_{20}P_4$  (free acid)

**MW** 1252.84 (free acid)

**Purity** > 95% HPLC

**Quantity** 0.5µmol = ~0.6mg

**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)