

# 3'-O-Piv-dATP

In pppApA a 5'-ATP unit is connected with a 5'-AMP unit via a 3' → 5' linkage to form a linear dinucleotide. pppApA is considered to be a linear dinucleotide intermediate in the enzymatic production of the bacterial signalling nucleotide c-diAMP (Cat. No. C 088) and of cyclic oligoadenylylates such as c-tetraAMP (Cat. No. C 335).

*BLG-T069-05 (5 x 1 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.*

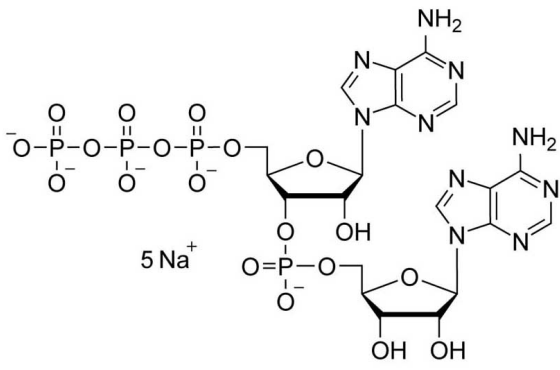
## Ordering Information

[Order Online »](#)

BLG-T069-01	1µmol
BLG-T069-05	5x1µmol

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

Alternative Name 5'-Triphosphoadenylyl-(3'→5')-adenosine . sodium salt

CAS 17730-98-2

Couple Type Modified nucleotides

Formula  $C_{20}H_{28}N_{10}O_{19}P_4$

Formulation 10 mM aqueous solution

MW 836.4

Purity > 95% HPLC

Quantity 1µmol is equivalent to ~0.8 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)