

# 3'-O-Ac-dATP

## Ribose-modified dATP analogue

3'-O-Ac-dATP is an analogue of dATP with blocked 3'-hydroxy function, which can be used as a transient, pH-sensitive chain terminator during DNA polymerization. The acetyl group of 3'-O-Ac-dATP is sensitive towards alkaline conditions and can be easily split off to allow further subsequent polymerization steps.

*BLG-A276-25 (5 x 5 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.*

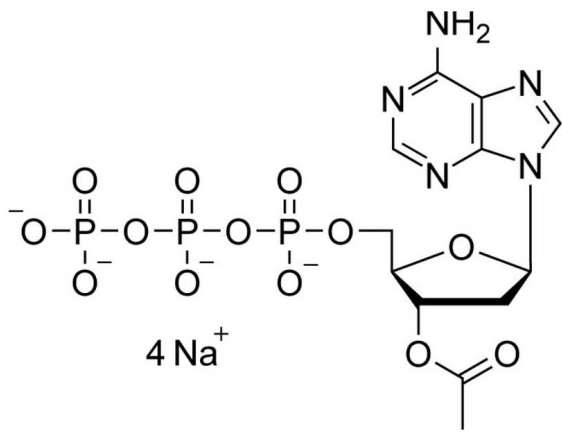
### Ordering Information

[Order Online »](#)

BLG-A276-05	5µmol
BLG-A276-25	5x5µmol

### 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** 3'-O-Acetyl-2'-deoxyadenosine-5'-O-triphosphate . sodium salt

**CAS** 28120-63-0

**Formula**  $C_{12}H_{18}N_5O_{13}P_3 \cdot 4 Na$

**Formulation** 10 mM aqueous solution

**MW** 533.2 (free acid)

**Purity** > 95% HPLC



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)