

# Ap<sub>4</sub> / Up<sub>4</sub>A

Dinucleoside polyphosphate containing both a purine nucleobase as well as a pyrimidine nucleobase moiety. Considered to be a novel endothelium-derived vasoconstrictive factor.

*BLG-U008-05 (5 x 1 µmol pack size) is not sold in the U.S. or Canada.  
Please [contact us](#) for available options.*

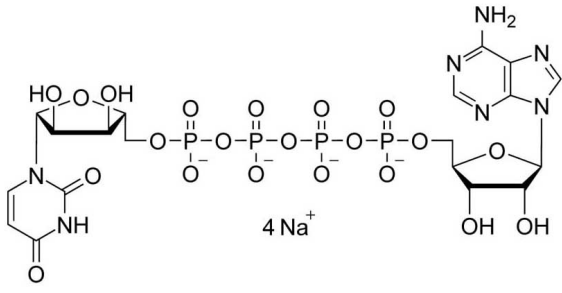
## Ordering Information

[Order Online »](#)

BLG-U008-05	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** P1-(5'-Adenosyl)-P4-(5'-uridy)-tetraphosphate . tetrasodium salt

**CAS** 10527-48-7

**Couple Type** Modified nucleotides

**Formula**  $C_{19}H_{23}N_7O_{21}P_4 \cdot 4Na$

**MW** 813.4 (free acid)

**Purity** > 95% HPLC

**Quantity** 1µmol≈0.8mg

**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)