

# ApCpp (10 mg)

## Ordering Information

[Order Online »](#)

JBS-NU-421-10	10mg
---------------	------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** AMPCPP, Adenosine-5'-[( $\alpha,\beta$ )-methylene]triphosphate, Sodium salt,  $\alpha,\beta$ -Methyleneadenosine-5'-triphosphate

**Appearance** white to off-white solid

**CAS** 7292-42-4 (acid), 1343364-54-4 (Trisodium salt)

**Couple Type** Modified nucleotides

**Formula**  $C_{11}H_{18}N_5O_{12}P_3$  (free acid)

**MW** 505.21 (free acid)

**Purity**  $\geq 98\%$  (HPLC)

**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-fr@enzolifesciences.com](mailto:info-fr@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)