

# ATP $\gamma$ S (5 mg)

Agonistic ligand, mainly for nucleoside receptor A1. Nucleosidephosphates stabilized against hydrolytic degradation can directly bind to nucleoside receptors.

## Ordering Information

[Order Online »](#)

JBS-NU-406-5	5mg
--------------	-----

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage -20°C

Shipping Dry Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** Adenosine-5'-(γ-thio)-triphosphate, Tetralithium salt, Adenosine-5'-(3-thio)-triphosphate, Adenosine-5'-(3-thiotriphosphate)

**Appearance** white to off-white solid

**Application Notes** Modulation of intracellular signalling  
Signalling of purinergic receptors  
Regulation of cation channels  
Modulation of cytokine secretion  
Substrate for kinases  
5' end labeling with T4 Polynucleotide Kinase (T4 PNK) of DNA and RNA

**CAS** 93839-89-5

**Couple Type** Modified nucleotides

**Formula**  $C_{10}H_{16}N_5O_{12}P_3S$  (free acid)

**MW** 523.24 (free acid)

**Purity** ≥ 90 % (HPLC), contains < 10 % ADP

**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)