

ATP γ S (25 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-406-25	25mg
---------------	------

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

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com