

ATP α S, L pack

Ordering Information

[Order Online »](#)

JBS-NU-408L	5x25 μ l
-------------	--------------

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 Adenosine-5'-(α -thio)-triphosphate, Sodium salt; (1 : 1 Mixture of Rp and Sp isomers)

Appearance colorless to slightly yellow solution

Application Notes
Crystal structure in complex with adenylyl cyclase
Activation of beta-cell adenylyl cyclase
Potent agonist for P2Y11-receptor
Full agonist for rP2Y1-receptor

CAS 29220-54-0

Couple Type Modified nucleotides

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

Formulation solution in water

MW 523.24 (free acid)

Purity $\geq 95\%$ (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-
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)