

# EDA-ATP $\gamma$ S – ATTO-550

## Ordering Information

[Order Online »](#)

JBS-NU-1609-550	50 $\mu$ l
-----------------	------------

## 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** 2'/3'-O-(2-Aminoethyl-carbamoyl)-adenosine-5'-(γ-thio)-triphosphate, labeled with ATTO 550, Triethylammonium salt

**Appearance** red-violet solution

**Couple Type** Modified nucleotides

**Emission Maximum** 576 nm

**Excitation Maximum** 554 nm

**Formula**  $C_{50}H_{63}N_{10}O_{16}P_3S$  (free acid)

**Formulation** solution in water

**MW** 1185.09 (free acid)

**Purity** ≥ 90 % (HPLC), contains approx. 5 % EDA-ADP-dye

**Technical Info / Product** For the Original Manufacturer's data sheet please [click here](#).

### Notes



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)