

EDA-GTP γ S – ATTO-565

Ordering Information

[Order Online »](#)

| | |
|-----------------|------------|
| JBS-NU-1610-565 | 50 μ 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)-guanosine-5'-(γ-thio)-triphosphate, labeled with ATTO 565, Triethylammonium salt

Couple Type Modified nucleotides

Emission Maximum 590 nm

Excitation Maximum 564 nm

Formula $C_{44}H_{50}N_9O_{18}P_3S$ (free acid)

Formulation solution in water

MW 1117.91 (free acid)

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

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