

# GTP $\gamma$ S (10 mg)

## Ordering Information

[Order Online »](#)

JBS-NU-412-10	10mg
---------------	------

## 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** Guanosine-5'-(γ-thio)-triphosphate, Tetralithium salt, Guanosine-5'-(3-thio)-triphosphate, Guanosine-5'-(3-thiotriphosphate)

**Appearance** white to off-white solid

**Application Notes** Functional assay for GPCRs  
Binding assay

**CAS** 94825-44-2

**Couple Type** Modified nucleotides

**Formula**  $C_{10}H_{16}N_5O_{13}P_3S$  (free acid)

**MW** 539.25 (free acid)

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

**Quantity** Unit definition: 1 unit = 1 µl of a 10 mM solution.

**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)