

GTP γ S (20 mg)

Ordering Information

[Order Online »](#)

JBS-NU-412-20	20mg
---------------	------

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

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