

# Rp-GTP- $\alpha$ -S

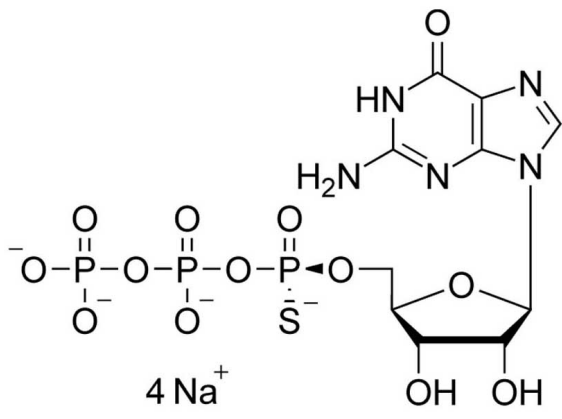
Inhibitor or activator of GTP binding proteins with often increased metabolic stability. Inhibitor of photoreceptor guanylate cyclase. Useful for characterization of GTP-responsive receptors and determination of their stereospecificity.

*BLG-G014-25 (5 x 5  $\mu$ mol pack size) is not sold in the U.S. or Canada.  
Please [contact us](#) for available options.*

## Ordering Information [Order Online »](#)

|                    |               |
|--------------------|---------------|
| <b>BLG-G014-05</b> | 5 $\mu$ mol   |
| <b>BLG-G014-25</b> | 5x5 $\mu$ mol |

## Manuals, SDS & CofA [View Online »](#)



## Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** Guanosine-5'-O-(1-thiotriphosphate) . tetrasodium salt, Rp-isomer

**Appearance** 10 mM aqueous solution

**CAS** 81570-50-5

**Couple Type** Modified nucleotides

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

**MW** 539.3 (free acid)

**Purity** > 95% HPLC

**Quantity** 5µmol≈2.7mg

**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)