

# TNP- $\beta$ -L-GTP

## Fluorescent analogue of $\beta$ -L-GTP

Fluorescent analogue of  $\beta$ -L-GTP, the mirror image version of naturally occurring  $\beta$ -D-GTP, where the guanine nucleobase is connected to  $\beta$ -L-ribose. TNP fluorophore with  $\lambda_{exc}$  408 nm and  $\lambda_{em}$  552 nm. It is sensitive to indicators of local environment such as polarity and viscosity, and frequently exhibits a spectral shift and fluorescence enhancement upon binding to a protein. Potential inhibitor of various adenylyl cyclase isoforms and soluble guanylyl cyclase (sGC).

*BLG-G035-05 (5 x 1  $\mu$ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.*

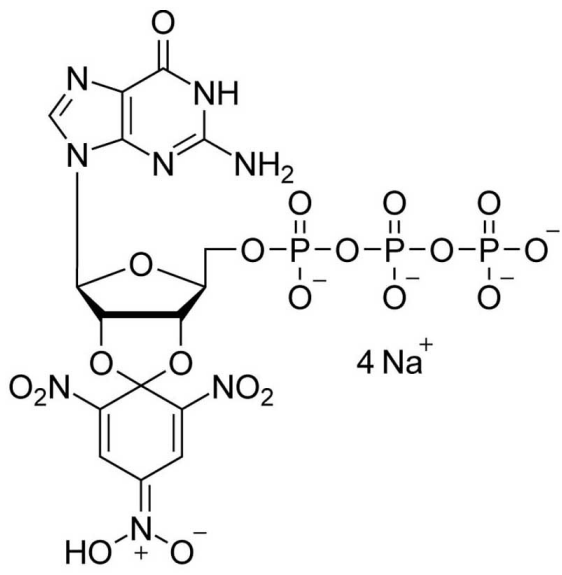
### Ordering Information

[Order Online »](#)

BLG-G035-01	1 $\mu$ mol
BLG-G035-05	5x1 $\mu$ 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** Guanine-2',3'-O-(2, 4, 6-trinitrophenyl)-β-L-ribofuranosyl-5'-O-triphosphate . sodium salt

**Couple Type** Modified nucleotides

**Formula**  $C_{16}H_{17}N_8O_{20}P_3$

**MW** 734.3

**Purity** > 95% HPLC

**Quantity** 1μmol is equivalent to ~0.7 mg

**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)