

# 3'-NH<sub>2</sub>-GTP

Ribose-modified GTP analogue

3'-NH<sub>2</sub>-GTP is an analogue of guanosine-5'-triphosphate, where the ribose 3' hydroxy group has been replaced by an amino group, which allows for conjugation with amine-reactive ligands to yield e.g. fluorescent or immobilized products.

*BLG-A250-25 (5 x 5 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.*

## Ordering Information

[Order Online »](#)

BLG-A250-05

5µmol

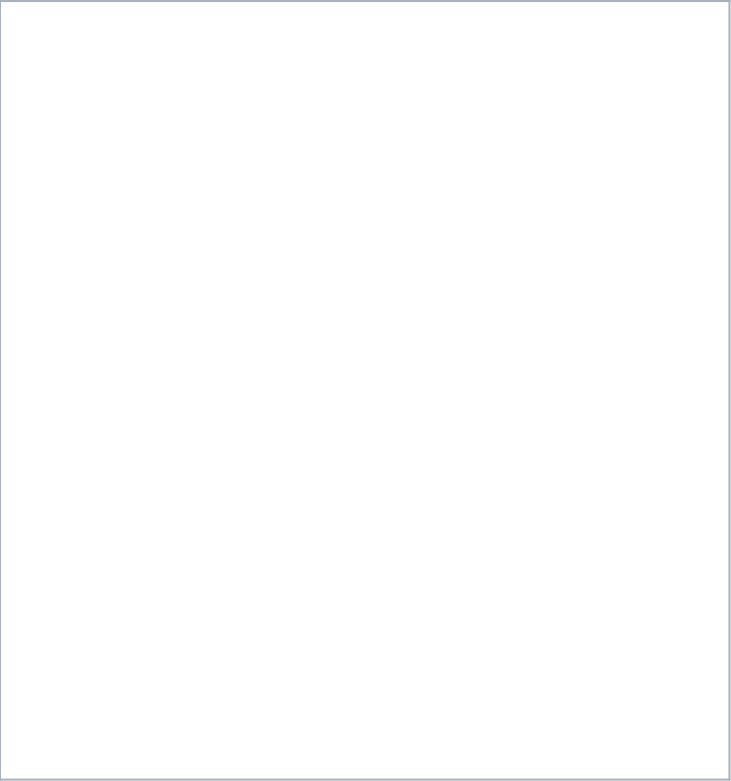
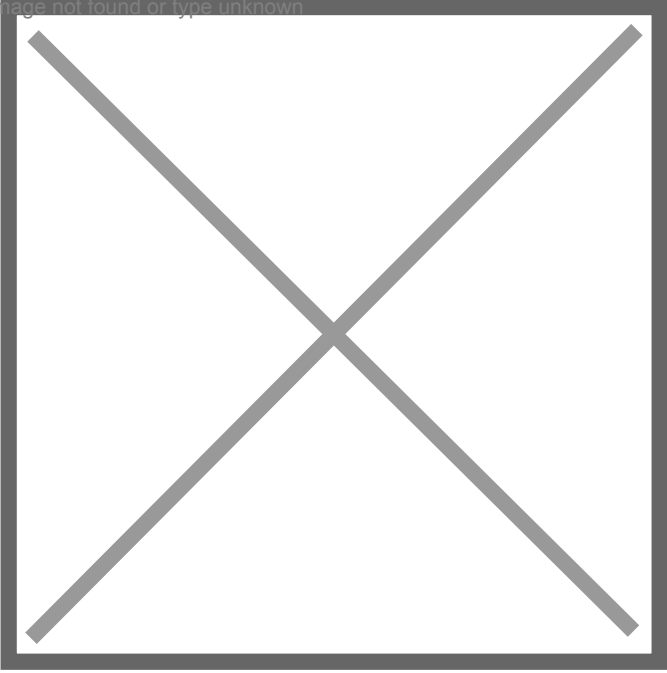
BLG-A250-25

5x5 $\mu$ mol

**Manuals, SDS & CofA**

[View Online »](#)

Image not found or type unknown



# Handling & Storage

Long Term Storage -80°C

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** 3'-Amino-3'-deoxyguanosine-5'-O-triphosphate . sodium salt

**CAS** 344402-44-4

**Couple Type** Modified nucleotides

**Formula**  $C_{10}H_{17}N_6O_{13}P_3$  (free acid)

**Formulation** 10 mM aqueous solution

**MW** 522.2 (free acid)

**Purity** > 95% HPLC

**Quantity** 5µmol is equivalent to ~2.6 mg

**Technical Info / Product Notes** For the Original Manufacturer's data sheet please [click here](#).

Image not found or type not known  
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)