

3'-NH₂-GTP

Ribose-modified GTP analogue

3'-NH₂-GTP is an analogue of guanosine-5'-triphosphate, where the ribose 3' hydroxy group has been replaced by an amino group, which allows for conjugation with amine-reactive ligands to yield e.g. fluorescent or immobilized products.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.*

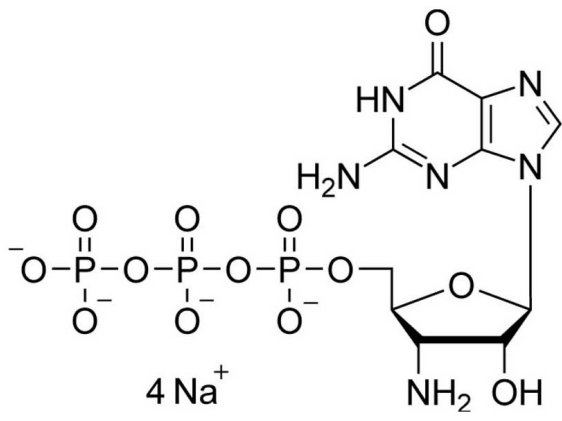
Ordering Information

[Order Online »](#)

BLG-A250-05	5µmol
BLG-A250-25	5x5µmol

Manuals, SDS & CofA

[View Online »](#)



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](#).



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