

β -L-GTP

Ribose-modified GTP analogue

β -L-GTP is the mirror image version of naturally occurring β -D-GTP, where the guanine nucleobase is connected to β -L-ribose. Nucleoside triphosphate precursor for the formation of highly nuclease-resistant L-RNA. L-RNA is used and explored for biological applications like aptamers and drug nanocarriers. Useful for research into GTP-dependent receptor proteins.

BLG-G030-25 (5 x 5 μ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

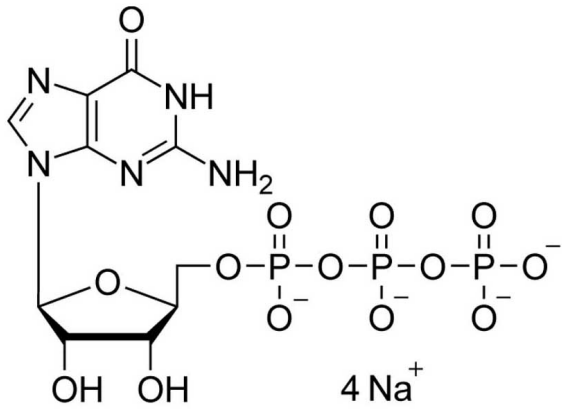
Ordering Information

[Order Online »](#)

BLG-G030-05	5 μ mol
BLG-G030-25	5x5 μ 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-β-L-ribofuranosyl-5'-O-triphosphate

CAS 784151-35-5

Couple Type Modified nucleotides

Formula $C_{10}H_{16}N_5O_{14}P_3$ (free acid)

MW 523.2 (free acid)

Purity > 95% HPLC

Quantity 5μmol are 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