

2'-O-Bn-3'-MANT-GTP

Ribose-modified fluorescent GTP analogue

2'-O-Bn-3'-MANT-GTP is a fluorescent (λ_{exc} 350 nm; λ_{em} 442 nm) analogue of GTP where the ribose hydroxy groups have been modified by an ether bond to a benzyl group and an ester bond to a fluorescent MANT group, respectively. Significantly improved stability against esterases compared to other 2'/3'-esters of the ribose like MANT-GTP (Cat. No. M 032). Potential inhibitor of various adenylyl cyclase isoforms and soluble guanylyl cyclase (sGC).

BLG-B187-05 (5 x 1 μ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

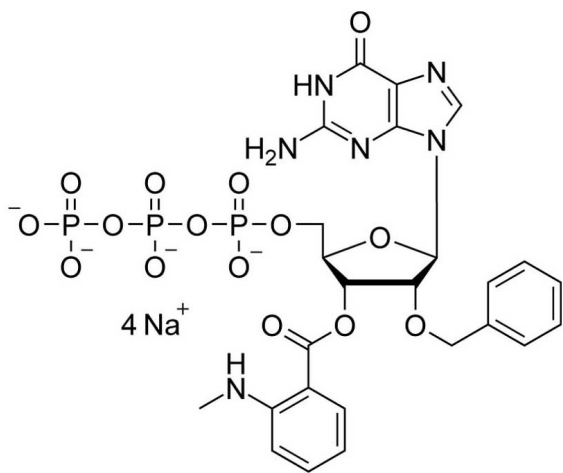
Ordering Information

[Order Online »](#)

BLG-B187-01	1 μ mol
BLG-B187-05	5x1 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 2'-O-Benzyl-3'-O-(N'-methylantraniloyl)guanosine-5'-O-triphosphate . sodium salt

Couple Type Modified nucleotides

Formula $C_{25}H_{29}N_6O_{15}P_3$ (free acid)

Formulation 1 mM aqueous solution

MW 746.5 (free acid)

Purity > 95% HPLC

Quantity 1 µmol is equivalent to ~0.7 mg



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)