

2'-O-Bn-GTP

Ribose-modified GTP analogue

2'-O-Bn-GTP is an analogue of GTP where the ribose 2'-hydroxy group has been modified by an ether bond to a benzyl group. Significantly improved stability against esterases compared to other 2'/3'-esters of the ribose like MANT-GTP (Cat. M 032). Potential inhibitor of various adenylyl cyclase isoforms and soluble guanylyl cyclase (sGC).

BLG-B183-05 (5 x 1 μ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

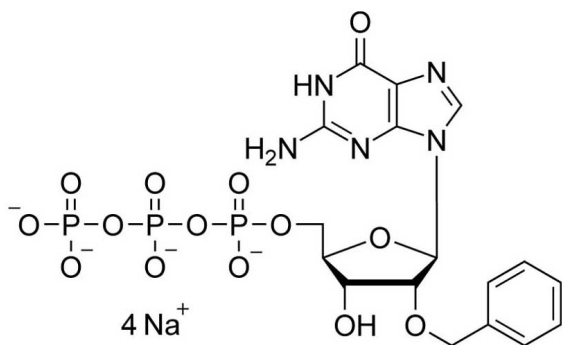
Ordering Information

[Order Online »](#)

BLG-B183-01	1 μ mol
BLG-B183-05	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-Benzylguanosine-5'-O-triphosphate . sodium salt

CAS 1610778-74-9

Couple Type Modified nucleotides

Formula $C_{17}H_{22}N_5O_{14}P_3$ (free acid)

Formulation 10 mM aqueous solution

MW 613.3 (free acid)

Purity > 95% HPLC

Quantity 1 μ mol is equivalent to ~0.6 mg



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