

8-Br-ITP

Analogue of ATP and GTP, respectively

8-Br-ITP is an analogue of inosine- 5'-O- triphosphate where the hydrogen atom in position 8 of the purine nuclepbase has been replaced by bromine.

*BLG-B157-05 (5 x 1 μ mol pack size) is not sold in the U.S. or Canada.
Please [contact us](#) for available options.*

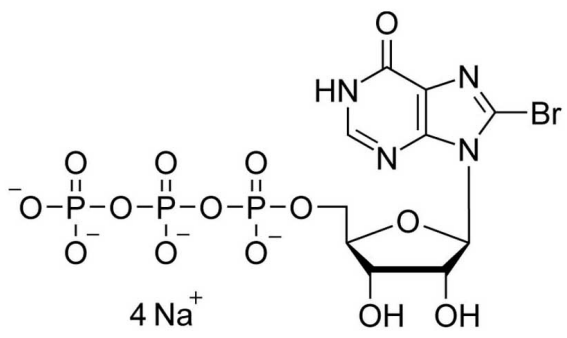
Ordering Information

[Order Online »](#)

BLG-B157-01	1 μ mol
BLG-B157-05	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 8-Bromoinosine-5'-O-triphosphate . sodium salt

CAS 81481-67-6

Couple Type Modified nucleotides

Formula $C_{10}H_{14}BrN_4O_{14}P_3$

Formulation 10 mM aqueous solution

MW 587.1

Purity > 95% HPLC

Quantity 1µmol is equivalent to ~0.6 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