

ANT-GTP

ANT-GTP is an analogue of guanosine-5'-triphosphate, where either the ribose 2' or the 3' hydroxyl group has been modified with the relatively small fluorescent anthraniloyl group (λ_{exc} 330 nm, λ_{em} ~428nm). The ANT fluorophore has a certain sensitivity for its environment and can change its spectral properties upon binding. Inhibitor of various adenylyl cyclase isoforms.

BLG-A214-25 (5 x 5 μ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

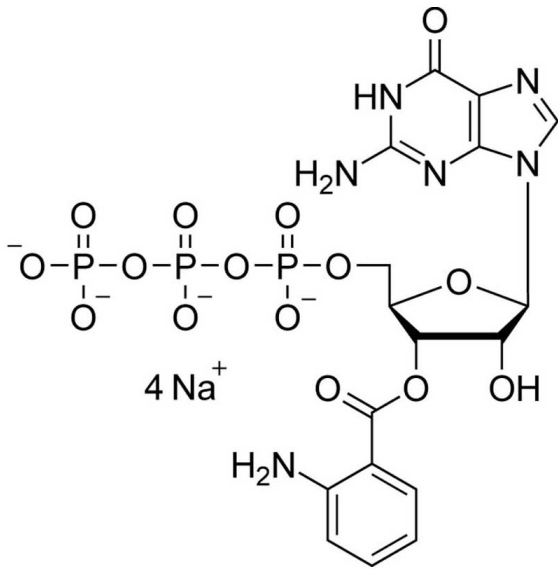
Ordering Information

[Order Online »](#)

BLG-A214-05	5 μ mol
BLG-A214-25	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 2'-/3'-O-Anthraniloylguanosine-5'-O-triphosphate . sodium salt

Couple Type Modified nucleotides

Formula $C_{17}H_{21}N_6O_{15}P_3$

MW 642.31

Purity > 95% HPLC

Quantity 5µmol are equivalent to ~3.2 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