

CI-ANT-ITP

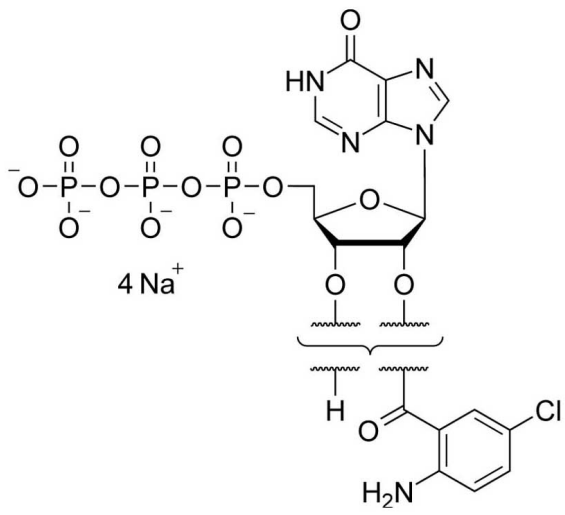
Analogue of inosine-5'-triphosphate with only moderate intrinsic fluorescence (λ_{exc} 350 nm; λ_{em} ~450 nm), increasing considerably within hydrophobic environment. Inhibitor of mammalian adenylate cyclases.

BLG-C132-25 (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-C132-05	5 μ mol
BLG-C132-25	5x5 μ mol

Manuals, SDS & CofA [View Online »](#)



Handling & Storage

Handling Protect from light.

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 2'-/3'-O-(5-Chloroanthraniloyl)inosine-5'-O-triphosphate . tetrasodium salt

Appearance 10 mM aqueous solution

CAS 1350521-48-0

Couple Type Modified nucleotides

Formula $C_{17}H_{19}ClN_5O_{15}P_3$ (free acid)

MW 661.7

Purity > 95% HPLC



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