

2-MeS-TNP-ITP

Fluorescent ITP analogue

Fluorescent analogue of ATP and GTP, respectively with λ_{exc} 408 nm and λ_{em} 552 nm, which is supposed to substitute for ATP and GTP in the interaction with various enzymes and proteins. Potential agonist of P2Y purinoceptors and supposed to be inhibitor of soluble guanylate cyclase. It is sensitive to indicators of local environment such as polarity and viscosity, and frequently exhibits a spectral shift and fluorescence enhancement upon binding to a protein.

BLG-M124-05 (5 x 1 μ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

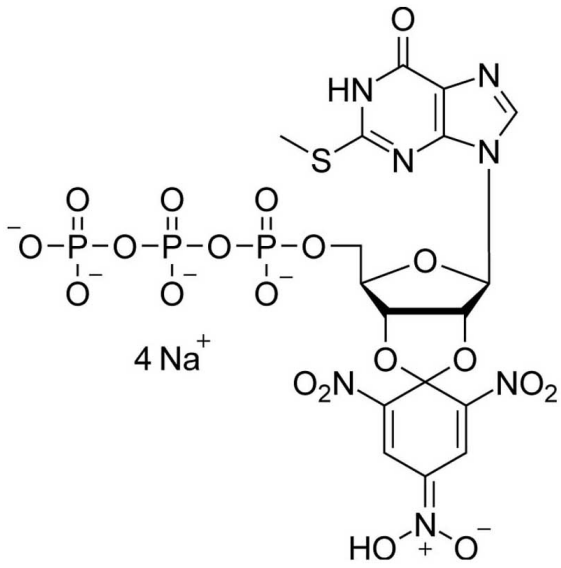
Ordering Information

[Order Online »](#)

BLG-M124-01	1 μ mol
BLG-M124-05	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-Methylthio-2',3'-O-(2,4,6-trinitrophenyl)inosine-5'-O-triphosphate . sodium salt

Couple Type Modified nucleotides

Formula $C_{17}H_{18}N_7O_{20}P_3S$

MW 765.4

Purity > 95% HPLC

Quantity 1µmol is equivalent to ~0.8 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