

2-MeS-ITP

Potential P2Y purinoceptor agonist

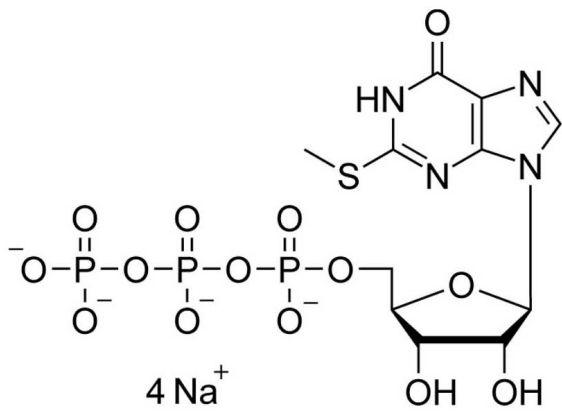
2-MeS-ITP is an analogue and potential metabolite of 2-MeS-ATP (Cat. No. M 021) where the 6-amino group has been replaced by an oxygen atom. Potential agonist of P2Y purinoceptors and supposed to be an inhibitor of soluble guanylate cyclase.

BLG-M114-05 (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-M114-01	1 μ mol
BLG-M114-05	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-Methylthioinosine-5'-O-triphosphate

Couple Type Modified nucleotides

Formula $C_{11}H_{17}N_4O_{14}P_3S$ (free acid)

Formulation 10 mM aqueous solution

MW 554.3 (free acid)

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