

# MANT-ITP- $\gamma$ -S

## Fluorescent and phosphorothioate-modified analogue of ITP

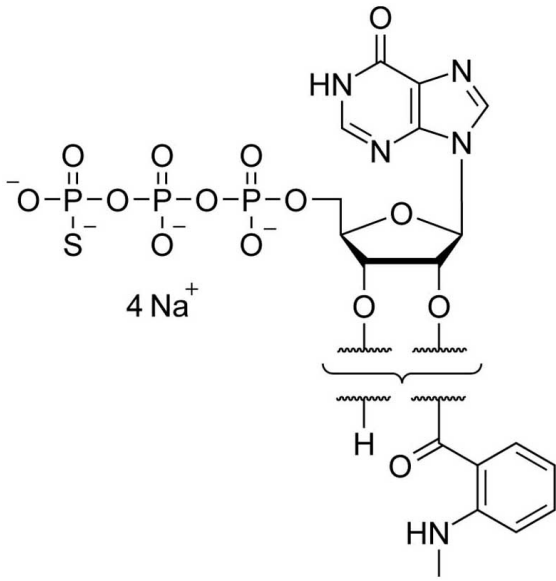
MANT-ITP- $\gamma$ -S is a fluorescent and phosphorothioate-modified analogue of the purine metabolism intermediate ITP with increased metabolic stability. The MANT fluorophore has a certain sensitivity for its environment and can change its spectral properties upon binding.

*BLG-M173-05 (5 x 1  $\mu$ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.*

### Ordering Information [Order Online »](#)

BLG-M173-01	1 $\mu$ mol
BLG-M173-05	5x1 $\mu$ 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-(N-Methylanthraniloyl)inosine-5'-O-(3-thiotriphosphate) . sodium salt

**Couple Type** Modified nucleotides

**Formula**  $C_{18}H_{22}N_5O_{14}P_3S$

**Formulation** 10 mM aqueous solution

**MW** 657.4

**Purity** > 95% HPLC for mixture

**Quantity** 1µmol is equivalent to ~0.7 mg



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)