

ITP- γ -S

Phosphorothioate-modified analogue of ITP

ITP- γ -S is a phosphorothioate-modified analogue of ITP with increased metabolic stability which can be useful for modification with SH-reactive reporters or for connection to structures with SH-groups via a disulfide bond. The parent compound ITP is an intermediate in the purine metabolism pathway. Gille et al. (2004) reported ITP- γ -S to be a mixed G-protein activator and adenylyl cyclase (AC) inhibitor.

*BLG-I046-05 (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-I046-01	1 μ mol
BLG-I046-05	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 Inosine-5'-O-(3-thiotriphosphate) . sodium salt

CAS 77065-87-3

Couple Type Modified nucleotides

Formula $C_{10}H_{15}N_4O_{13}P_3S$

Formulation 10 mM aqueous solution

MW 524.2

Purity > 95% HPLC

Quantity 0.1 µmol are equivalent to ~0.5 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