

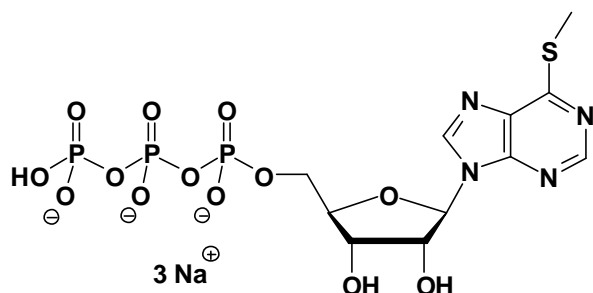
6-Methylthio-ITP

6-Methylthioinosine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-1133S	150 Units
NU-1133L	750 Units

Selected References:

Lavi *et al.* (1985) A rapid and sensitive high-performance liquid chromatographic assay for 6-mercaptopurine metabolites in red blood cells. *Anal. Biochem.* **144**(2):514.



Cat. No.: NU-1133

Molecular Formula: $C_{11}H_{14}N_4O_{13}P_3S$ (Anion)

Molecular Weight: 535.23 (Anion)

Purity: > 95%, clear aqueous solution, pH 7.5

Spectroscopic properties: λ_{max} 224/292 nm;
 ϵ 11500/19000

Storage conditions:

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at < -20°C. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

For research use only!

1 unit = 1 μ l of a 10 mM solution