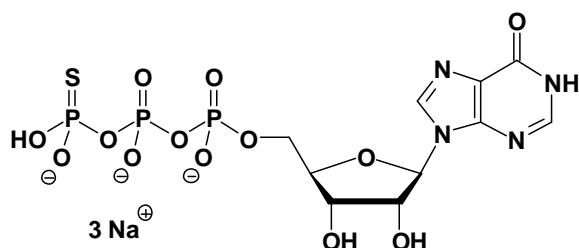


ITP γ S

Inosine-5'-(γ -thio)-triphosphate, Sodium salt

Cat. No.	Amount
NU-413S	20 Units
NU-413L	100 Units



Cat. No.: NU-413

Molecular Formula: $C_{10}H_{12}N_4O_{13}P_3S$ (Anion)

Molecular Weight: 521.20 (Anion)

Purity: > 90%, clear aqueous solution, pH 7.5

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 6 months after date of delivery.

For research use only!

1 unit = 1 μ l of a 10 mM solution