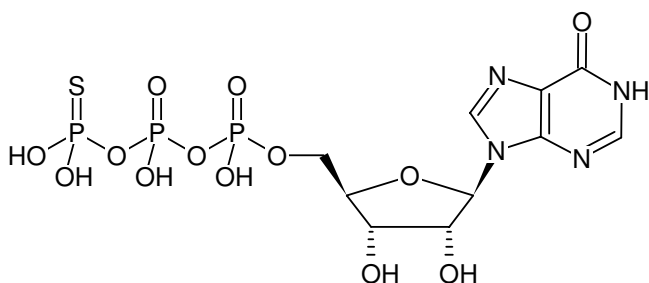


**ITPyS**

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

Cat. No.	Amount
NU-413S	20 µl (10 mM)
NU-413L	5 x 20 µl (10 mM)



Structural formula of ITPyS

For general laboratory use.**Shipping:** shipped on dry ice**Storage Conditions:** store at -20 °C**Shelf Life:** 6 months after date of delivery**Molecular Formula:** C₁₀H₁₅N₄O₁₃P₃S (free acid)**Molecular Weight:** 524.23 g/mol (free acid)**Exact Mass:** 523.96 g/mol (free acid)**CAS#:** 77065-87-3**Purity:** ≥ 90 % (HPLC)**Form:** solution in water**Color:** colorless to slightly yellow**Concentration:** 10 mM - 11 mM**pH:** 7.5 ±0.5**Spectroscopic Properties:** λ_{max} 249 nm, ε 12.2 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.0)