

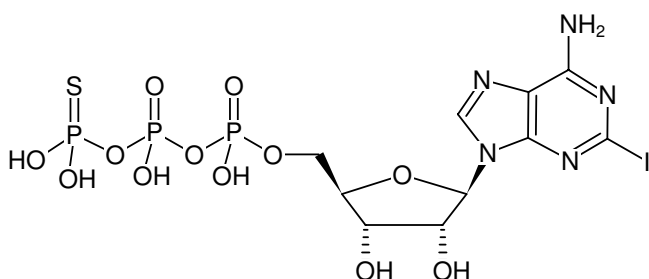
**2-Iodo-ATPyS**

(2I-ATPyS)

2-Iodo-adenosine-5'-(γ-thio)-triphosphate, Sodium salt

Cat. No.	Amount
NU-905-1	1 mg
NU-905-5	5 mg

Please note: For reasons of stability, please make sure that the pH value of a solution of this product never drops below 7.0. This can be achieved by dissolving the nucleotide in a buffer of your choice (50 - 100 mM, pH 7 - 10). Dissolve and adjust concentration photometrically.



Structural formula of 2-Iodo-ATPyS

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:** C₁₀H₁₅IN₅O₁₂P₃S (free acid)**Molecular Weight:** 649.14 g/mol (free acid)**Exact Mass:** 648.87 g/mol (free acid)**Purity:** ≥ 85 % (HPLC)**Form:** solid**Color:** white to off-white**Spectroscopic Properties:** λ_{max} 267 nm, ε 13.2 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)