

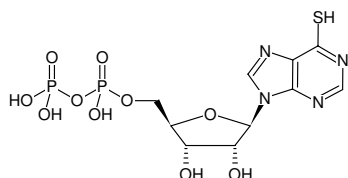


6-Mercaptopurine-riboside-5'-diphosphate

6-Thio-inosine-5'-diphosphate

6-Mercaptopurine-riboside-5'-diphosphate, Sodium salt

Cat. No.	Amount
NU-1210S	100 µl (10 mM)
NU-1210L	5 x 100 µl (10 mM)



Structural formula of 6-Mercaptopurine-riboside-5'-diphosphate

For research use only!

Shipping: shipped on gel packs

Storage Conditions: store at -20 °C

Short term exposure (up to 1 week cumulative) to ambient temperature possible.

Shelf Life: 12 months after date of delivery

Molecular Formula: C₁₀H₁₄N₄O₁₀P₂S (free acid)

Molecular Weight: 444.25 g/mol (free acid)

Exact Mass: 443.99 g/mol (free acid)

CAS#: 805-63-0

Purity: ≥ 95 % (HPLC)

Form: solution in water

Color: colorless to slightly yellow

Concentration: 10 mM - 11 mM

pH: 7.5 ±0.5

Spectroscopic Properties: λ_{max} 322 nm, ε 27.3 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)