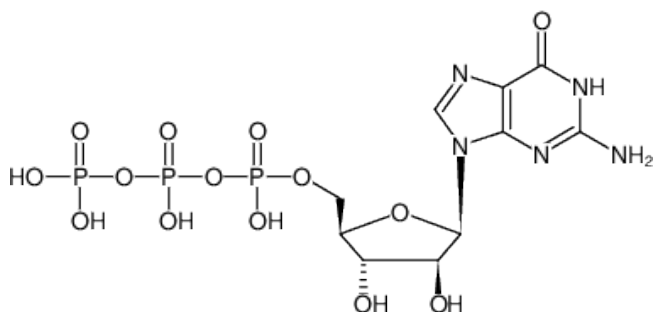



ara-Guanosine-5'-triphosphate (ara-GTP)

Guanosine-arabinofuranoside-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-1217S	50 µl (10 mM)
NU-1217L	5 x 50 µl (10 mM)



Structural formula of ara-Guanosine-5'-triphosphate (ara-GTP)

For general laboratory use.
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: C₁₀H₁₆N₅O₁₄P₃ (free acid)

Molecular Weight: 523.18 g/mol (free acid)

Exact Mass: 522.99 g/mol (free acid)

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} 252 nm, ε 13.6 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)