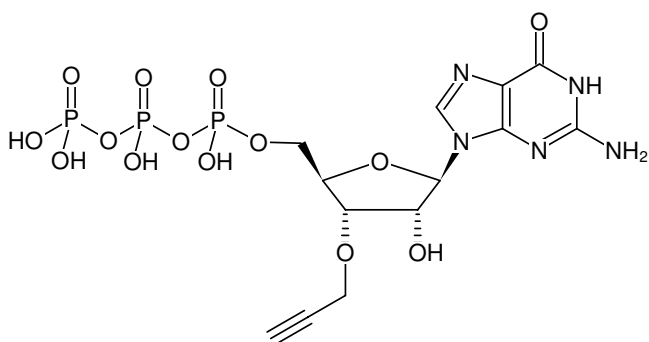




3'-(O-Propargyl)-GTP

3'-(O-Propargyl)-guanosine-5'-triphosphate, Triethylammonium salt

Cat. No.	Amount
NU-946	1 μ mol



Structural formula of 3'-(O-Propargyl)-GTP

For general laboratory use.

Shipping: shipped on gel packs

Storage Conditions: store at -20 °C

Additional Storage Conditions: Short term exposure (up to 1 week cumulative) to ambient temperature possible.

Shelf Life: 12 months after date of delivery

Molecular Formula: C₁₃H₁₈N₅O₁₄P₃ (free acid)

Molecular Weight: 561.23 g/mol (free acid)

Exact Mass: 561.01 g/mol (free acid)

CAS#: 1610778-63-6 (free acid)

Purity: \geq 95 % (HPLC)

Form: solid

Color: colorless to white

Spectroscopic Properties: λ_{max} 252 nm, ϵ 13.7 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)