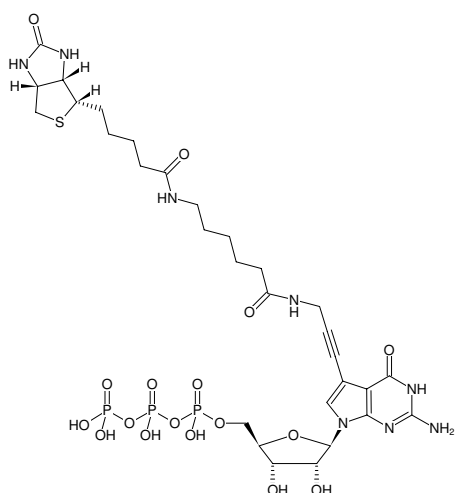


**Biotin-11-GTP**

Triethylammonium salt

Cat. No.	Amount
NU-971-BIOX-S	10 µl (1 mM)
NU-971-BIOX-L	5 x 10 µl (1 mM)



Structural formula of Biotin-11-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:** C₃₀H₄₅N₈O₁₇P₃S (free acid)**Molecular Weight:** 914.71 g/mol (free acid)**Exact Mass:** 914.18 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** filtered solution (30 kDa) in 10 mM Tris-HCl**Color:** colorless to slightly yellow**Concentration:** 1.0 mM - 1.1 mM**pH:** 7.5 ±0.5**Spectroscopic Properties:** λ_{max} 272 nm, ε 11.9 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)