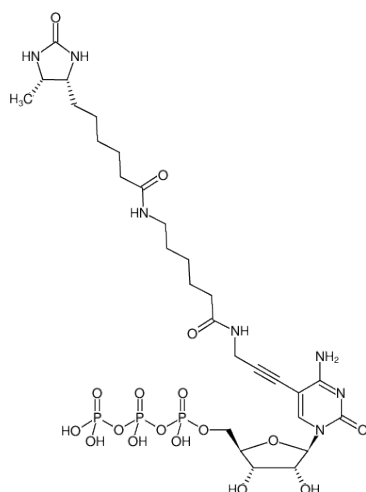




Desthiobiotin-11-CTP

5-Propargylamino-cytidine-5'-triphosphate, labeled with Desthiobiotin, Triethylammonium salt

| Cat. No. | Amount |
|-------------------|---------------|
| NU-831-Desthiobio | 30 µl (10 mM) |



Structural formula of Desthiobiotin-11-CTP

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: C₂₈H₄₆N₇O₁₇P₃ (free acid)

Molecular Weight: 845.63 g/mol (free acid)

Exact Mass: 845.22 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: filtered solution (30 kDa) in 10 mM Tris-HCl

Color: colorless to slightly yellow

Concentration: 10 mM – 11 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{max} 294 nm, ε 9.3 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)