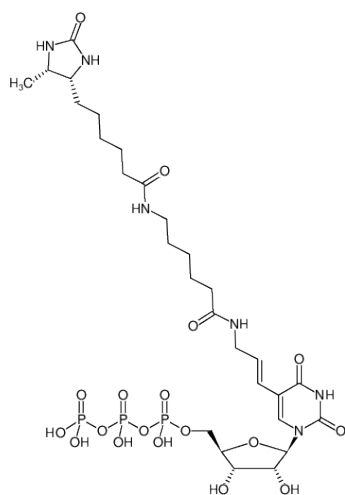


**Desthiobiotin-11-UTP**

5-(3-Aminoallyl)-uridine-5'-triphosphate, labeled with Desthiobiotin, Triethylammonium salt

Cat. No.	Amount
NU-821-Desthiobio	30 µl (10 mM)



Structural formula of Desthiobiotin-11-UTP

**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:** C<sub>28</sub>H<sub>47</sub>N<sub>6</sub>O<sub>18</sub>P<sub>3</sub> (free acid)**Molecular Weight:** 848.63 g/mol (free acid)**Exact Mass:** 848.22 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<sub>max</sub> 289 nm, ε 7.1 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)