

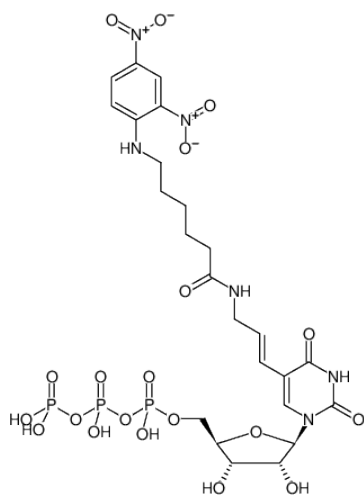


### Aminoallyl-UTP-DNP

Aminoallyl-UTP - Dinitrophenol, DNP-11-UTP

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

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



Structural formula of Aminoallyl-UTP-DNP

#### 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>24</sub>H<sub>33</sub>N<sub>6</sub>O<sub>20</sub>P<sub>3</sub>

**Molecular Weight:** 818.47 g/mol

**Exact Mass:** 818.10 g/mol

**Purity:** ≥ 95 % (HPLC)

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

**Color:** yellow

**Concentration:** 10 mM - 11 mM

**pH:** 7.5 ±0.5

**Spectroscopic Properties:** λ<sub>max</sub> 361 nm, ε 17.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)