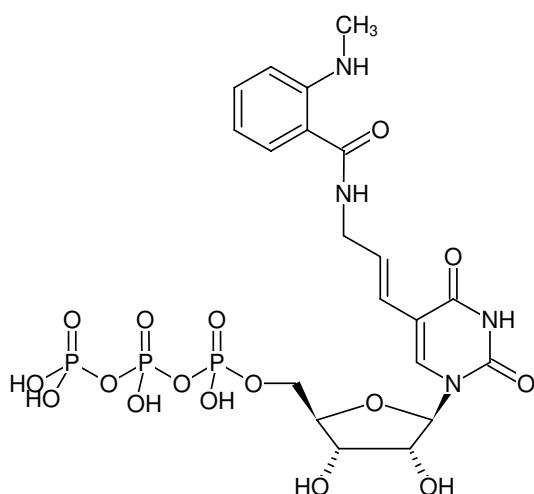


**Aminoallyl-UTP-MANT**

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

| Cat. No.   | Amount        |
|------------|---------------|
| NU-821-MNT | 100 µl (5 mM) |



Structural formula of Aminoallyl-UTP-MANT

**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>20</sub>H<sub>27</sub>N<sub>4</sub>O<sub>16</sub>P<sub>3</sub> (free acid)**Molecular Weight:** 672.37 g/mol (free acid)**Exact Mass:** 672.06 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** solution in water**Color:** colorless to slightly yellow**Concentration:** 5.0 mM - 5.5 mM**pH:** 7.5 ±0.5**Spectroscopic Properties:** λ<sub>exc</sub> 335 nm, λ<sub>em</sub> 440 nm, ε 2.7 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)