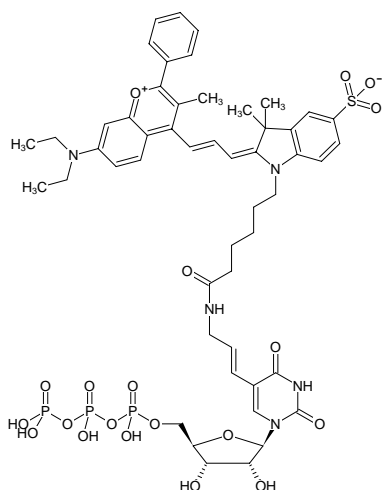




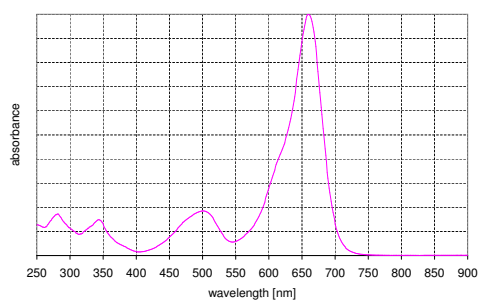
## Aminoallyl-UTP - DYQ-660

5-(3-Aminoallyl)-uridine-5'-triphosphate, labeled with DYQ 660, Triethylammonium salt

Cat. No.	Amount
NU-821-660Q	40 µl (1 mM)



Structural formula of Aminoallyl-UTP - DYQ-660



absorption spectrum of DYQ 660

### For research use only!

**Shipping:** shipped on blue ice

**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>51</sub>H<sub>62</sub>N<sub>5</sub>O<sub>20</sub>P<sub>3</sub>S (free acid)

**Molecular Weight:** 1190.05 g/mol (free acid)

**Purity:** ≥ 95 % (HPLC)

**Form:** clear aqueous solution

**Concentration:** 1 mM - 1.1 mM

**pH:** 7.5 ± 0.5

**Spectroscopic Properties:** λ<sub>exc</sub> 660 nm; ε 109.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)

λ<sub>exc</sub> 660 nm; ε 140.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Ethanol)