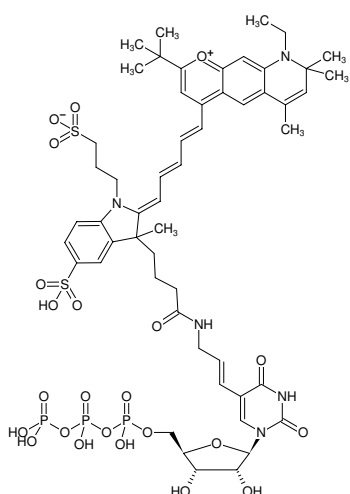




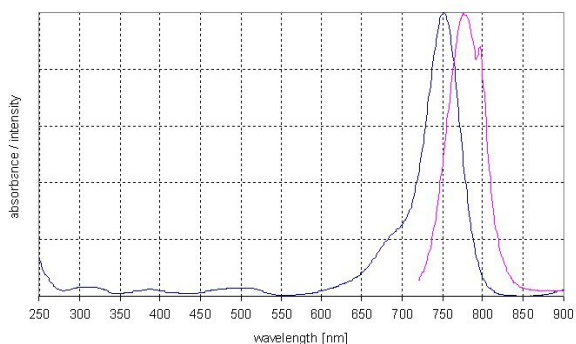
### Aminoallyl-UTP-DY-751

5-(3-Aminoallyl)-uridine-5'-triphosphate, labeled with DY 751, Triethylammonium salt

Cat. No.	Amount
NU-821-751	40 µl (1 mM)



Structural formula of Aminoallyl-UTP-DY-751



excitation and emission spectrum of DY 751

**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>54</sub>H<sub>70</sub>N<sub>5</sub>O<sub>23</sub>P<sub>3</sub>S<sub>2</sub> (free acid)

**Molecular Weight:** 1313.31 g/mol (free acid)

**Purity:** ≥ 95 % (HPLC)

**Form:** solution in water

**Concentration:** 1.0 mM - 1.1 mM

**pH:** 7.5 ± 0.5

**Spectroscopic Properties:** λ<sub>exc</sub> 751 nm, λ<sub>em</sub> 779 nm, ε 270.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Ethanol)