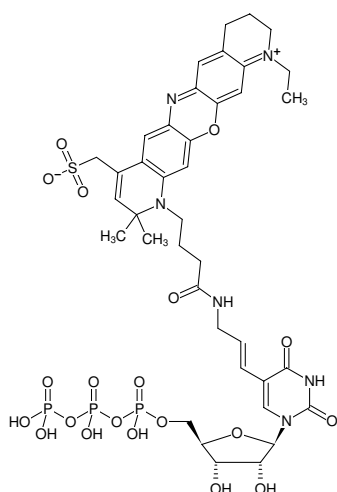




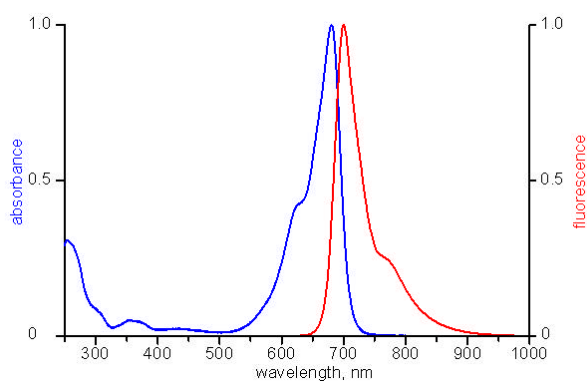
Aminoallyl-UTP-ATTO-680

5-(3-Aminoallyl)-uridine-5'-triphosphate, labeled with ATTO 680, Triethylammonium salt

Cat. No.	Amount
NU-821-680	10 µl (1 mM)



Structural formula of Aminoallyl-UTP-ATTO-680



excitation and emission spectrum of ATTO 680

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: C₃₉H₄₉N₆O₂₀P₃S (free acid)

Molecular Weight: 1046.82 g/mol (free acid)

Exact Mass: 1046.19 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_{exc} 681 nm, λ_{em} 698 nm, ε 125.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)

Applications:

Fluorescence-labeling of RNA^[1,2]

Analysis of RNA-protein interaction^[2]

Selected References:

[1] Besse *et al.* (2009) *Drosophila* PTB promotes formation of high-order RNP particles and represses *oskar* translation. *Genes Dev.* **23**:195.

[2] Köhn *et al.* (2010) Near-infrared (NIR) dye-labeled RNAs identify binding of ZBP1 to the noncoding Y3-RNA. *RNA* **16** (7):1420.