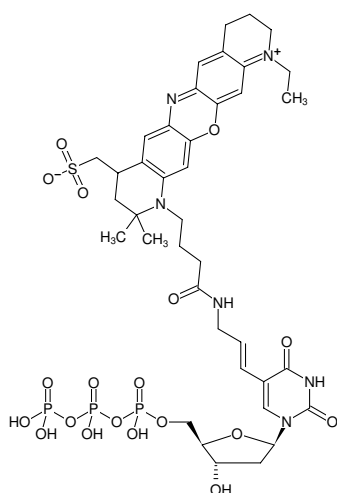




Aminoallyl-dUTP-ATTO-655

5-(3-Aminoallyl)-2'-deoxyuridine-5'-triphosphate, labeled with ATTO 655, Triethylammonium salt

Cat. No.	Amount
NU-803-655-S	10 µl (1 mM)
NU-803-655-L	5 x 10 µl (1 mM)



Structural formula of Aminoallyl-dUTP-ATTO-655

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: C₃₉H₅₁N₆O₁₉P₃S (free acid)

Molecular Weight: 1032.84 g/mol (free acid)

Exact Mass: 1032.21 g/mol (free acid)

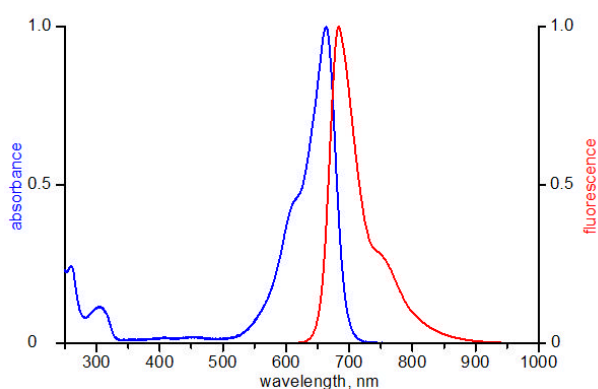
Purity: ≥ 95 % (HPLC)

Form: clear aqueous solution

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 663 nm, λ_{em} 680 nm, ε 125.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



excitation and emission spectrum of ATTO 655