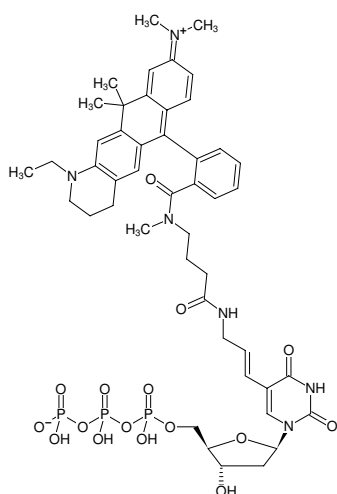




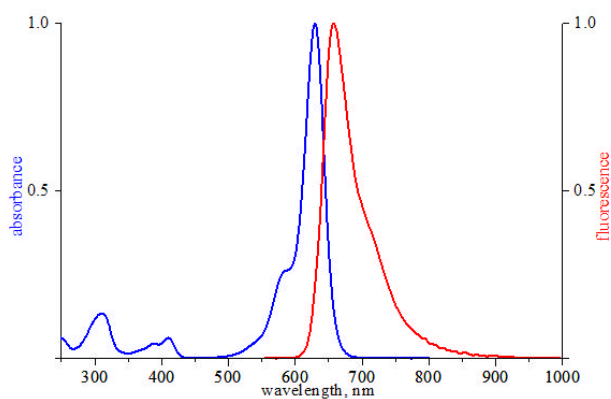
Aminoallyl-dUTP-ATTO-633

5-(3-Aminoallyl)-2'-deoxyuridine-5'-triphosphate, labeled with ATTO 633, Triethylammonium salt

Cat. No.	Amount
NU-803-633	40 µl (1 mM)



Structural formula of Aminoallyl-dUTP-ATTO-633



excitation and emission spectrum of ATTO 633

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: C₄₇H₅₉N₆O₁₆P₃ (free acid)

Molecular Weight: 1056.94 g/mol (free acid)

Exact Mass: 1056.32 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} 630 nm, λ_{em} 651 nm, ε 130.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)

Selected References:

Xiang *et al.* (2017) Correlative live and super-resolution imaging reveals the dynamic structure of replication domains *bioRxiv* doi:10.1101/189373.