



5-Propargylamino-ddUTP-ATTO-643

5-Propargylamino-2',3'-dideoxyuridine-5'-triphosphate, labeled with ATTO 643, Triethylammonium salt

Cat. No.	Amount
NU-1619-643	40 µl (1 mM)

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: C₁₂H₁₇N₃O₁₃P₃ - ATTO 643 (free acid)

Molecular Weight: 1323.26 g/mol (free acid)

Exact Mass: 1322.35 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

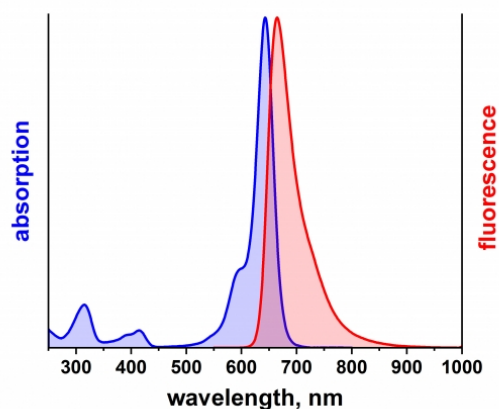
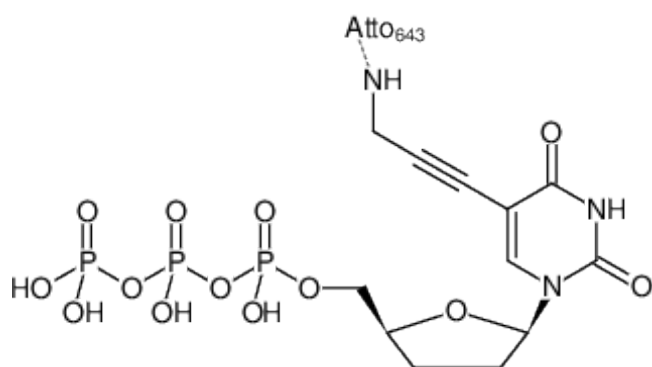
Form: solution in water

Color: blue

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 643 nm, λ_{em} 665 nm, ε 150.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



Excitation and Emission spectrum of ATTO 643