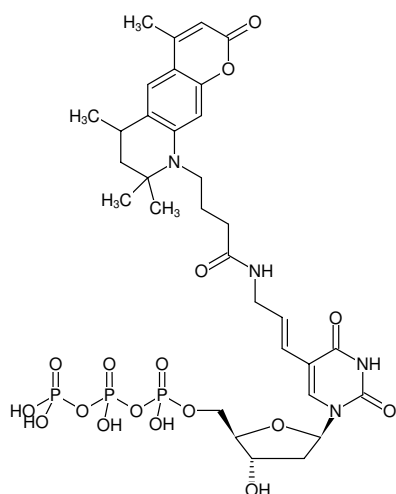




Aminoallyl-dUTP-ATTO-390

5-(3-Aminoallyl)-2'-deoxyuridine-5'-triphosphate, labeled with ATTO 390, Triethylammonium salt

Cat. No.	Amount
NU-803-390	80 µl (1 mM)



Structural formula of Aminoallyl-dUTP-ATTO-390

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: C₃₂H₄₃N₄O₁₇P₃ (free acid)

Molecular Weight: 848.63 g/mol (free acid)

Exact Mass: 848.18 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

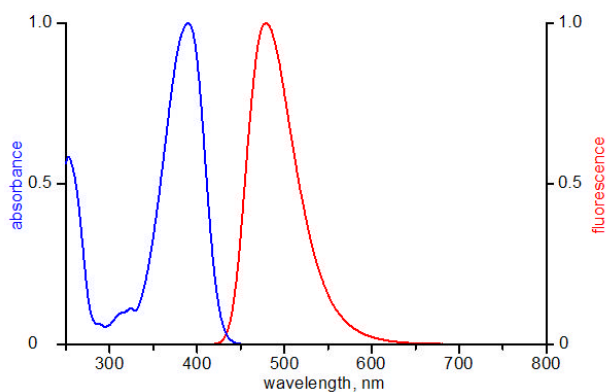
Form: solution in water

Color: slightly yellow

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 390 nm, λ_{em} 476 nm, ε 24.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



excitation and emission spectrum of ATTO 390