



Aminoallyl-dUTP-ATTO-612Q

5-(3-Aminoallyl)-2'-deoxyuridine-5'-triphosphate, labeled with ATTO 612Q, Triethylammonium salt

Cat. No.	Amount
NU-803-612Q	40 μ l (1 mM)

For general laboratory use.

Shipping: shipped on gel packs

Storage Conditions: store at $-20\text{ }^{\circ}\text{C}$

Short term exposure (up to 1 week cumulative) to ambient temperature possible.

Shelf Life: 12 months after date of delivery

Molecular Formula: $\text{C}_{12}\text{H}_{20}\text{N}_3\text{O}_{14}\text{P}_3$ - ATTO 612Q (free acid)

Molecular Weight: 1195.22 g/mol (free acid)

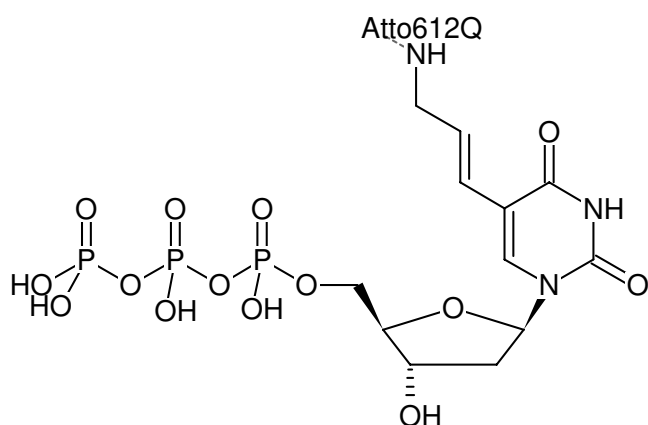
Purity: $\geq 95\%$ (HPLC)

Form: solution in water

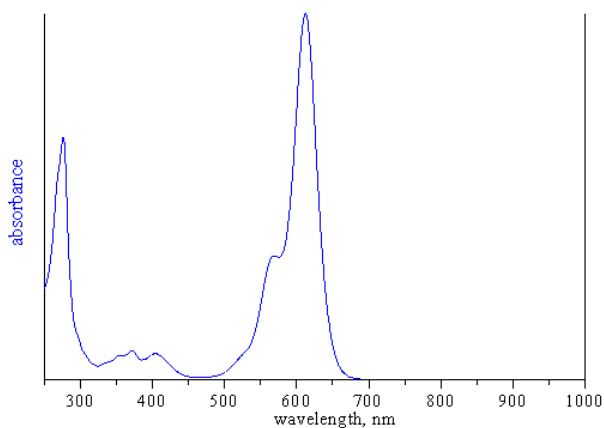
Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 615 nm, ϵ 115.0 L $\text{mmol}^{-1}\text{ cm}^{-1}$ (Tris-HCl pH 7.5)



Structural formula of Aminoallyl-dUTP-ATTO-612Q



absorption spectrum of ATTO 612Q