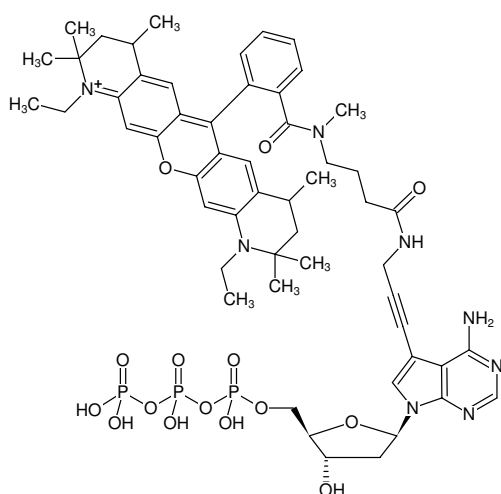




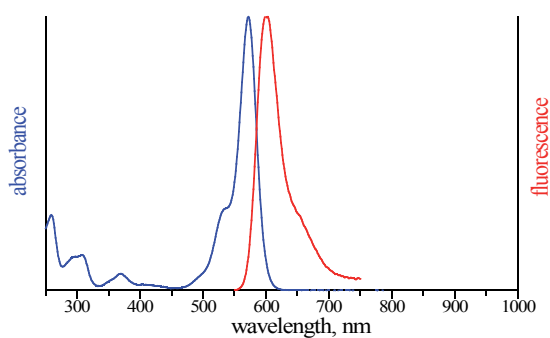
7-Propargylamino-7-deaza-ATTO-Rho12

7-Deaza-7-propargylamino-2'-deoxyadenosine-5'-triphosphate, labeled with ATTO Rho12, Triethylammonium salt

| Cat. No. | Amount |
|---------------|--------------|
| NU-1611-RHO12 | 40 µl (1 mM) |



Structural formula of 7-Propargylamino-7-deaza-dATP-ATTO-Rho12



excitation and emission spectrum of ATTO Rho12

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: C₅₅H₆₉N₈O₁₅P₃ (free acid)

Molecular Weight: 1175.12 g/mol (free acid)

Exact Mass: 1174.41 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: solution in water

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 576 nm, λ_{em} 601 nm, ε 120.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)