

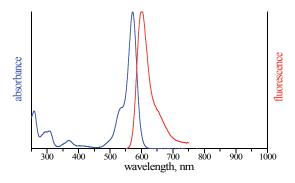


■ 7-Propargylamino-7-deaza-ddATP-ATTO-Rho12

7-Deaza-7-propargylamino-2',3'-dideoxyadenosine-5'-triphosphate, labeled with ATTO Rho12, Triethylammonium salt

Cat. No.	Amount
NU-1612-RHO12	80 μl (0,5 mM)

Structural formula of 7-Propargylamino-7-deaza-ddATP-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: $C_{55}H_{69}N_8O_{14}P_3$ (free acid) Molecular Weight: 1159.12 g/mol (free acid)

Exact Mass: 1158.41 g/mol (free acid)

Purity: ≥ 95 % (HPLC) **Form:** solution in water

Concentration: 0.50 mM - 0.55 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)