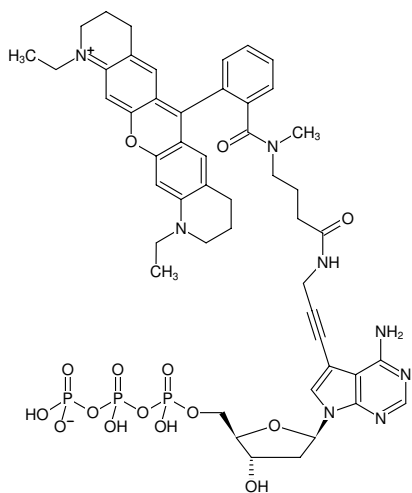




### 7-Propargylamino-7-deaza-dATP-ATTO-Rho11

7-Deaza-7-propargylamino-2'-deoxyadenosine-5'-triphosphate, labeled with ATTO Rho11, Triethylammonium salt

Cat. No.	Amount
NU-1611-RHO11	40 µl (1 mM)



Structural formula of 7-Propargylamino-7-deaza-dATP-ATTO-Rho11

**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<sub>49</sub>H<sub>57</sub>N<sub>8</sub>O<sub>15</sub>P<sub>3</sub> (free acid)

**Molecular Weight:** 1090.96 g/mol (free acid)

**Exact Mass:** 1090.32 g/mol (free acid)

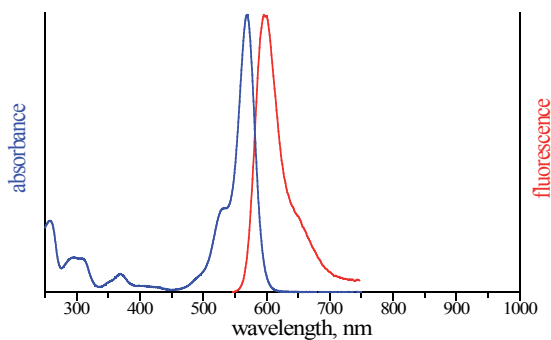
**Purity:** ≥ 95 % (HPLC)

**Form:** solution in water

**Concentration:** 1.0 mM - 1.1 mM

**pH:** 7.5 ± 0.5

**Spectroscopic Properties:** λ<sub>exc</sub> 572 nm, λ<sub>em</sub> 595 nm, ε 120.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)



excitation and emission spectrum of ATTO Rho11