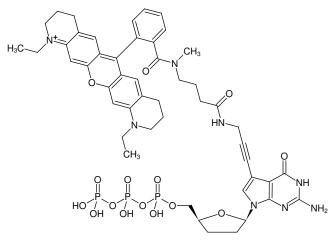




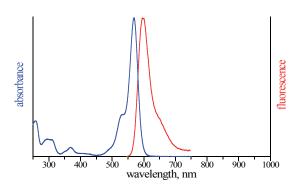
## 7-Propargylamino-7-deaza-ddGTP-ATTO-Rho11

7-Deaza-7-propargylamino-2',3'-dideoxyguanosine-5'-triphosphate, labeled with ATTO Rho11, Triethylammonium salt

Cat. No.	Amount
NU-1618-RHO11	80 μl (0,5 mM)



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



excitation and emission spectrum of 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:**  $C_{49}H_{57}N_8O_{15}P_3$  (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: 0.50 mM - 0.55 mM

pH: 7.5 ±0.5

Spectroscopic Properties:  $\lambda_{exc}$  572 nm,  $\lambda_{em}$  595 nm,

 $\epsilon$  120.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)