

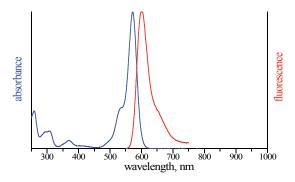


## 5-Propargylamino-ddUTP-ATTO-Rho12

5-Propargylamino-2',3'-dideoxyuridine-5'-triphosphate, labeled with ATTO Rho12, Triethylammonium salt

Cat. No.	Amount
NU-1619-RHO12	80 μl (0,5 mM)

 $Structural\ formula\ of\ 5-Propargy lamino-ddUTP-ATTO-Rho12$ 



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:**  $C_{53}H_{67}N_6O_{16}P_3$  (free acid) **Molecular Weight:** 1137.07 g/mol (free acid)

Exact Mass: 1136.38 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:  $\lambda_{exc}$  576 nm,  $\lambda_{em}$  601 nm,

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