



### 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)

**For research use only!**

**Shipping:** shipped on blue ice

**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>53</sub>H<sub>67</sub>N<sub>6</sub>O<sub>16</sub>P<sub>3</sub> (free acid)

**Molecular Weight:** 1136.38 g/mol (free acid)

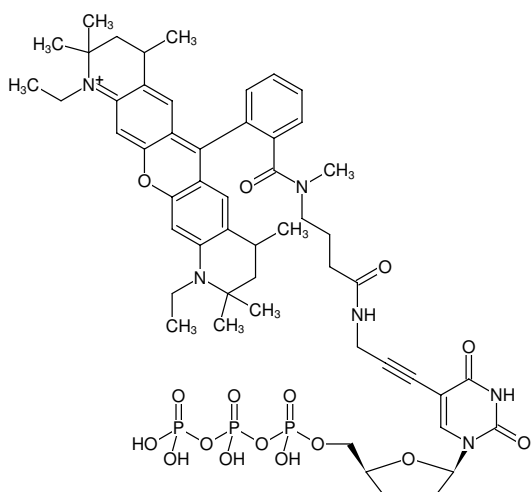
**Purity:** ≥ 95 % (HPLC)

**Form:** clear aqueous solution, pH 7.5 ± 0.5

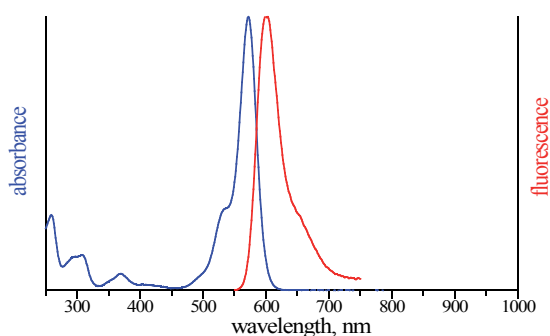
**Concentration:** 0.5 mM

**pH:** 7.5 ± 0.5

**Spectroscopic Properties:** λ<sub>exc</sub> 576 nm; λ<sub>em</sub> 601 nm; ε 120.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)



Structural formula of 5-Propargylamino-ddUTP - ATTO-Rho12



excitation and emission spectrum of ATTO Rho12