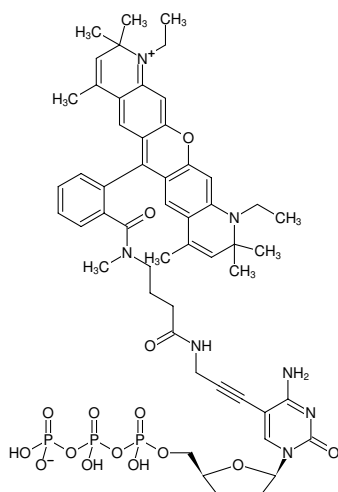




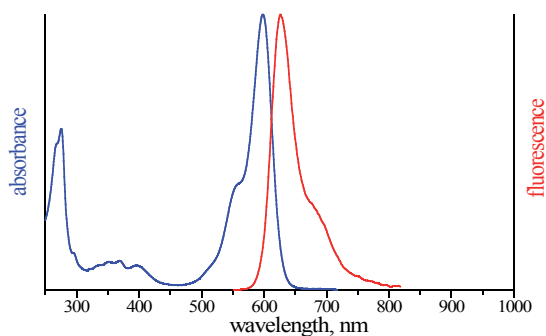
## 5-Propargylamino-ddCTP-ATTO-Rho13

5-Propargylamino-2',3'-dideoxycytidine-5'-triphosphate, labeled with ATTO Rho13, Triethylammonium salt

Cat. No.	Amount
NU-850-RHO13	80 $\mu$ l (0,5 mM)



Structural formula of 5-Propargylamino-ddCTP-ATTO-Rho13



excitation and emission spectrum of ATTO Rho13

For general laboratory use.

Please centrifuge briefly before opening (volume  $\leq$  2 ml).

**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>53</sub>H<sub>64</sub>N<sub>7</sub>O<sub>15</sub>P<sub>3</sub> (free acid)

**Molecular Weight:** 1132.05 g/mol (free acid)

**Exact Mass:** 1131.37 g/mol (free acid)

**Purity:**  $\geq$  95 % (HPLC)

**Form:** solution in water

**Color:** blue

**Concentration:** 0.50 mM - 0.55 mM

**pH:** 7.5  $\pm$  0.5

**Spectroscopic Properties:**  $\lambda_{exc}$  603 nm,  $\lambda_{em}$  627 nm,  $\epsilon$  120.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)