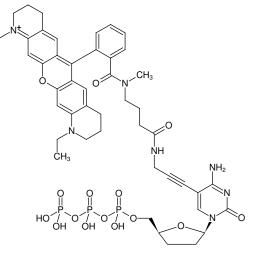




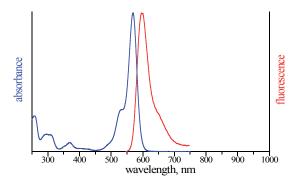
## 5-Propargylamino-ddCTP-ATTO-Rho11

5-Propargylamino-2',3'-dideoxycytidine-5'-triphosphate, labeled with ATTO Rho11, Triethylammonium salt

Cat. No.	Amount
NU-850-RHO11	80 μl (0,5 mM)



Structural formula of 5-Propargylamino-ddCTP-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<sub>47</sub>H<sub>56</sub>N<sub>7</sub>O<sub>15</sub>P<sub>3</sub> (free acid)

**Molecular Weight:** 1051.92 g/mol (free acid)

Exact Mass: 1051.30 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)

