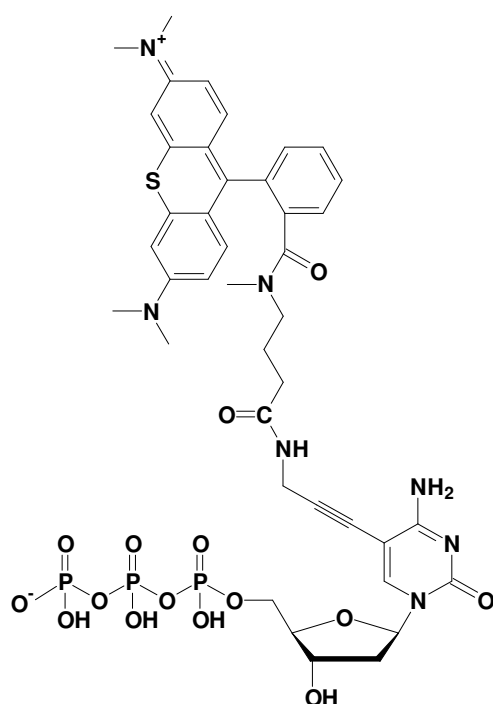


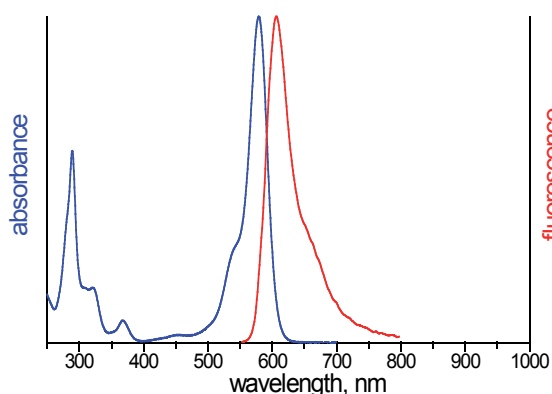
## 5-Propargylamino-dCTP - ATTO Thio12

5-Propargylamino-2'-deoxy-cytidine-5'-triphosphate, labeled with ATTO Thio12,  
 Triethylammonium salt

Cat. No.	Amount
NU-809-THIO12	40 $\mu$ l (1 mM)



structural formula of 5-Propargylamino-dCTP - ATTO Thio12



excitation and emission spectrum of ATTO Thio12

**Cat. No.:** NU-809-THIO12

**Molecular Formula:** C<sub>41</sub>H<sub>48</sub>N<sub>7</sub>O<sub>15</sub>P<sub>3</sub>S (free acid)

**Molecular Weight:** 1003,85 (free acid)

**Purity:** > 95%, clear aqueous solution, pH 7.5

**Spectroscopic properties:**

$\lambda_{exc}$  579 nm;  $\lambda_{em}$  609 nm;  $\epsilon$  1,1 x 10<sup>5</sup> cm<sup>-1</sup> M<sup>-1</sup>

**Storage conditions:**

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at < -20°C. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

**For research use only!**