



### 5-Propargylamino-ddUTP-ATTO-540Q

5-Propargylamino-2',3'-dideoxyuridine-5'-triphosphate, labeled with ATTO 540Q, Triethylammonium salt

Cat. No.	Amount
NU-1619-540Q	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>12</sub>H<sub>18</sub>N<sub>3</sub>O<sub>13</sub>P<sub>3</sub> - ATTO 540Q (free acid)

**Molecular Weight:** 1045.20 g/mol (free acid)

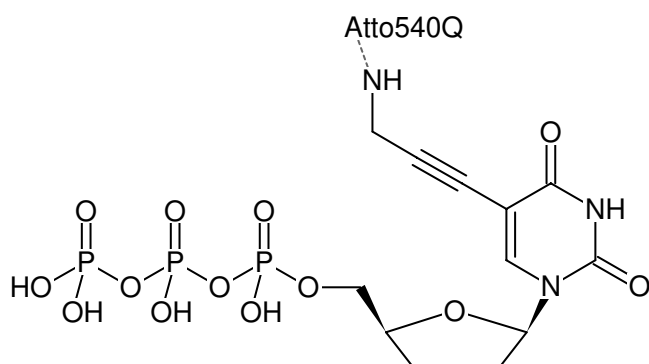
**Purity:** ≥ 95 % (HPLC)

**Form:** clear aqueous solution

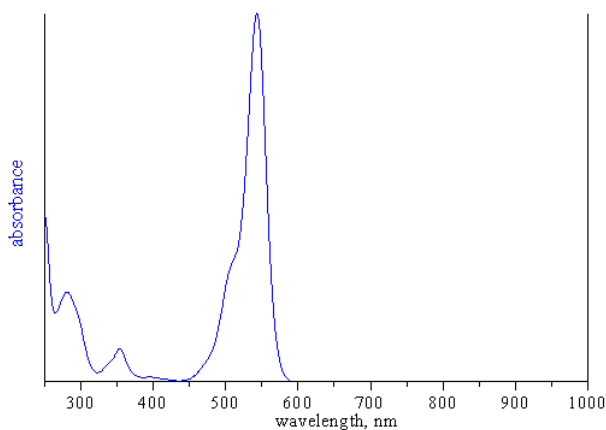
**Concentration:** 0.50 mM - 0.55 mM

**pH:** 7.5 ± 0.5

**Spectroscopic Properties:** λ<sub>exc</sub> 543 nm, ε 105.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)



Structural formula of 5-Propargylamino-ddUTP-ATTO-540Q



absorption spectrum of ATTO 540Q