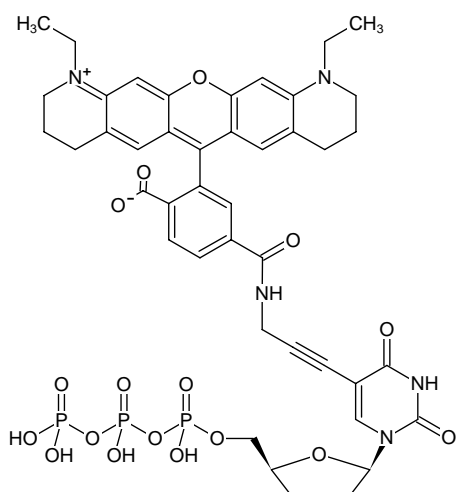




### 5-Propargylamino-ddUTP-ATTO-565

5-Propargylamino-2',3'-dideoxyuridine-5'-triphosphate, labeled with ATTO 565, Triethylammonium salt

| Cat. No.    | Amount       |
|-------------|--------------|
| NU-1619-565 | 80 µl (1 mM) |



Structural formula of 5-Propargylamino-ddUTP-ATTO-565

**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>43</sub>H<sub>46</sub>N<sub>5</sub>O<sub>17</sub>P<sub>3</sub> (free acid)

**Molecular Weight:** 997.78 g/mol (free acid)

**Exact Mass:** 997.21 g/mol (free acid)

**Purity:** ≥ 95 % (HPLC)

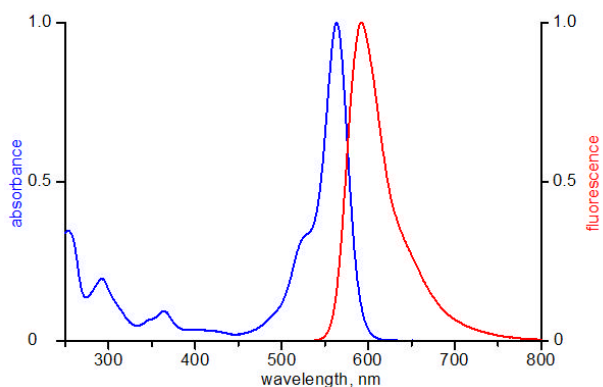
**Form:** solution in water

**Color:** red

**Concentration:** 1.0 mM - 1.1 mM

**pH:** 7.5 ± 0.5

**Spectroscopic Properties:** λ<sub>exc</sub> 564 nm, λ<sub>em</sub> 590 nm, ε 120.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)



excitation and emission spectrum of ATTO 565