



## 5-Propargylamino-ddUTP-DYQ-660

5-Propargylamino-2',3'-dideoxyuridine-5'-triphosphate, labeled with DYQ 660, Triethylammonium salt

Cat. No.	Amount
NU-1619-660Q	80 µl (0,5 mM)

**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>51</sub>H<sub>60</sub>N<sub>5</sub>O<sub>18</sub>P<sub>3</sub>S (free acid)

**Molecular Weight:** 1156.04 g/mol (free acid)

**Purity:** ≥ 95 % (HPLC)

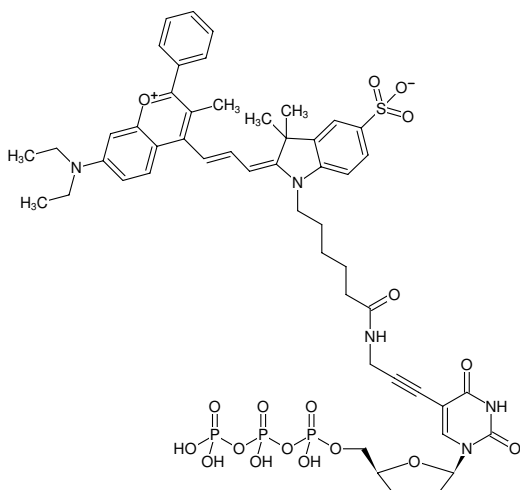
**Form:** solution in water

**Concentration:** 0.50 mM - 0.55 mM

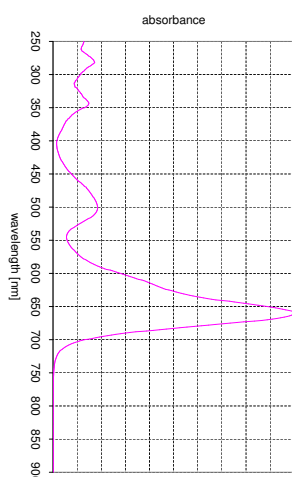
**pH:** 7.5 ± 0.5

**Spectroscopic Properties:** λ<sub>exc</sub> 660 nm, ε 109.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)

λ<sub>exc</sub> 660 nm, ε 140.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Ethanol)



Structural formula of 5-Propargylamino-ddUTP-DYQ-660



absorption spectrum of DYQ 660