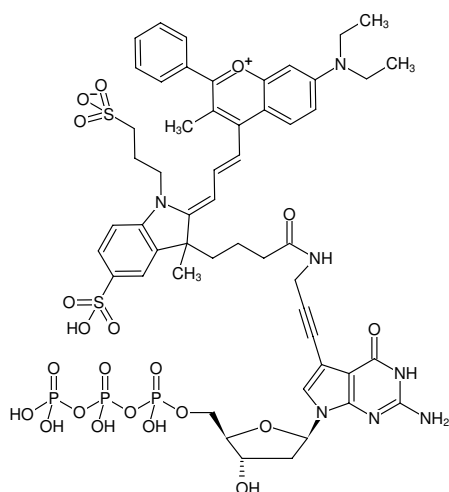




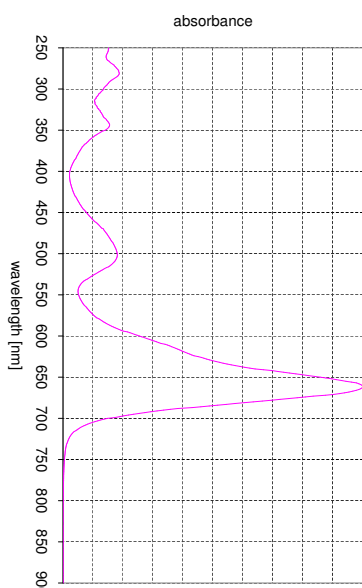
## 7-Propargylamino-7-deaza-dGTP-DYQ-661

7-Deaza-7-propargylamino-2'-deoxyguanosine-5'-triphosphate, labeled with DYQ 661, Triethylammonium salt

Cat. No.	Amount
NU-1615-661Q	40 µl (1 mM)



Structural formula of 7-Propargylamino-7-deaza-dGTP-DYQ-661



absorption spectrum of DYQ 661

**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>53</sub>H<sub>62</sub>N<sub>7</sub>O<sub>21</sub>P<sub>3</sub>S<sub>2</sub> (free acid)

**Molecular Weight:** 1290.15 g/mol (free acid)

**Exact Mass:** 1289.27 g/mol (free acid)

**Purity:** ≥ 95 % (HPLC)

**Form:** solution in water

**Concentration:** 1.0 mM - 1.1 mM

**pH:** 7.5 ± 0.5

**Spectroscopic Properties:** λ<sub>exc</sub> 662 nm, ε 104.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)

λ<sub>exc</sub> 662 nm, ε 140.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Ethanol)

**Please note:** DYQ661 is not a dark quencher, λ<sub>em</sub> 690 nm<sup>[1]</sup>

**Selected References:**

[1] Tan *et al.* (2010) Direct detection of Delta9-tetrahydrocannabinol in saliva using a novel homogeneous competitive immunoassay with fluorescence quenching. *Anal. Chim. Acta.* **658** (2):187.