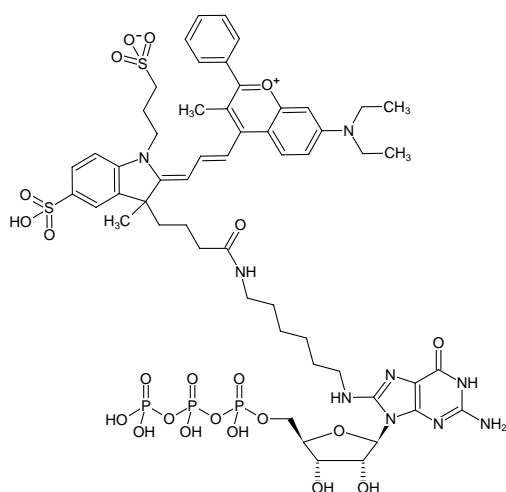




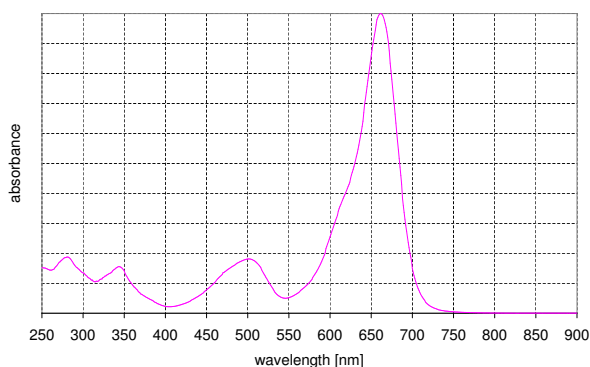
## 8-(6-Aminoethyl)-amino-GTP - DYQ-661

8-(6-Aminoethyl)-amino-guanosine-5'-triphosphate, labeled with DYQ 661, Triethylammonium salt

Cat. No.	Amount
NU-830-661Q	80 µl (1 mM)



Structural formula of 8-(6-Aminoethyl)-amino-GTP - DYQ-661



absorption spectrum of DYQ 661

### 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>55</sub>H<sub>71</sub>N<sub>9</sub>O<sub>22</sub>P<sub>3</sub>S<sub>2</sub> (Anion)

**Molecular Weight:** 1367.25 g/mol (Anion)

**Purity:** ≥ 95 % (HPLC)

**Form:** clear aqueous solution

**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.