



## $\gamma$ -(6-Aminoethyl)-ATP-DYQ-661

$\gamma$ -(6-Aminoethyl)-adenosine-5'-triphosphate, labeled with DYQ 661, Triethylammonium salt

Cat. No.	Amount
NU-833-661Q	80 $\mu$ l (1 mM)

For general laboratory use.

Please centrifuge briefly before opening (volume  $\leq$  2 ml).

**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>55</sub>H<sub>70</sub>N<sub>8</sub>O<sub>21</sub>P<sub>3</sub>S<sub>2</sub> (Anion)

**Molecular Weight:** 1336.24 g/mol (Anion)

**Purity:**  $\geq$  95 % (HPLC)

**Form:** solution in water

**Color:** blue

**Concentration:** 1.0 mM - 1.1 mM

**pH:** 7.5  $\pm$  0.5

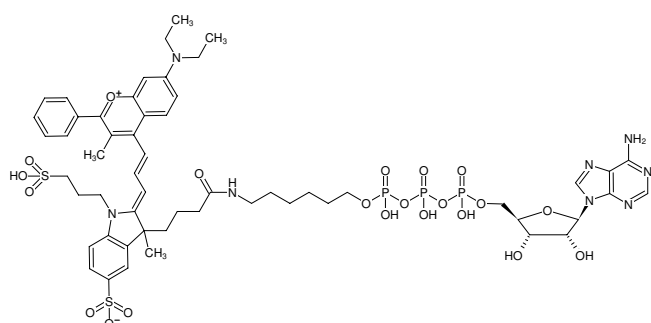
**Spectroscopic Properties:**  $\lambda_{exc}$  662 nm,  $\epsilon$  104.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)

$\lambda_{exc}$  662 nm,  $\epsilon$  140.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Ethanol)

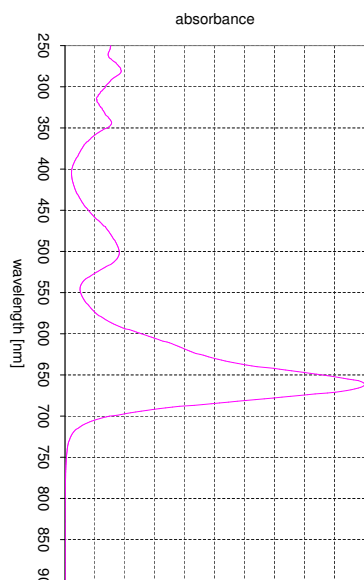
**Please note:** DYQ661 is not a dark quencher,  $\lambda_{em}$  690 nm<sup>[1]</sup>

**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.



Structural formula of  $\gamma$ -(6-Aminoethyl)-ATP-DYQ-661



absorption spectrum of DYQ 661