



EDA-ADP-DYQ-661

2'/3'-O-(2-Aminoethyl-carbamoyl)-Adenosine-5'-diphosphate, labeled with DYQ 661, Triethylammonium salt

Cat. No.	Amount
NU-802-661Q	40 µl (1 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: C₅₂H₆₂N₉O₁₉P₂S₂ (Anion)

Molecular Weight: 1243.18 g/mol (Anion)

Purity: ≥ 95 % (HPLC)

Form: solution in water

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

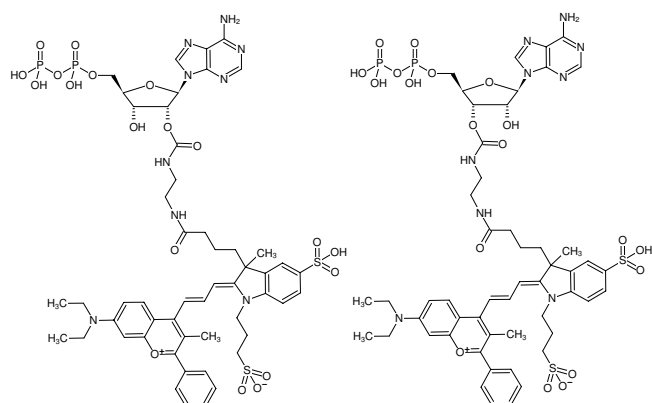
Spectroscopic Properties: λ_{exc} 662 nm, ε 104.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)

λ_{exc} 662 nm, ε 140.0 L mmol⁻¹ cm⁻¹ (Ethanol)

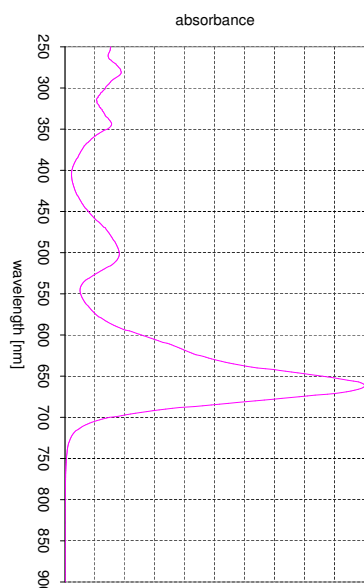
Please note: DYQ661 is not a dark quencher, λ_{em} 690 nm^[1]

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 EDA-ADP-DYQ-661



absorption spectrum of DYQ 661