



## EDA-ADP-DYQ-660

2'/3'-O-(2-Aminoethyl-carbamoyl)-adenosine-5'-diphosphate, labeled with DYQ 660, Triethylammonium salt

Cat. No.	Amount
NU-802-660Q	40 µl (1 mM)

For general laboratory use.

Please centrifuge briefly before opening (volume ≤ 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>52</sub>H<sub>63</sub>N<sub>9</sub>O<sub>16</sub>P<sub>2</sub>S (free acid)

**Molecular Weight:** 1164.13 g/mol (free acid)

**Purity:** ≥ 95 % (HPLC)

**Form:** solution in water

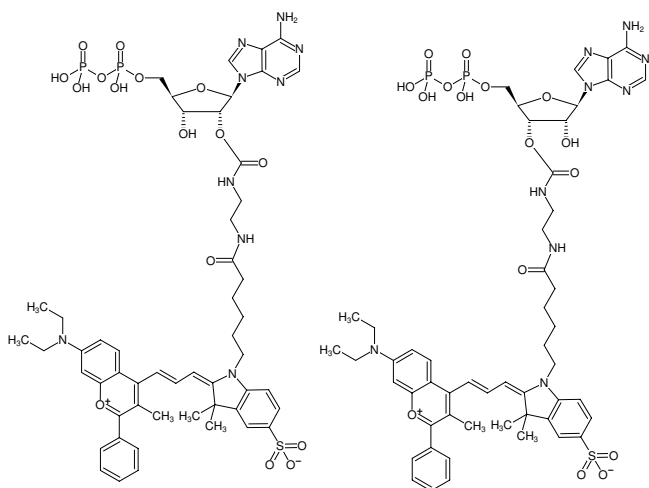
**Color:** blue

**Concentration:** 1.0 mM - 1.1 mM

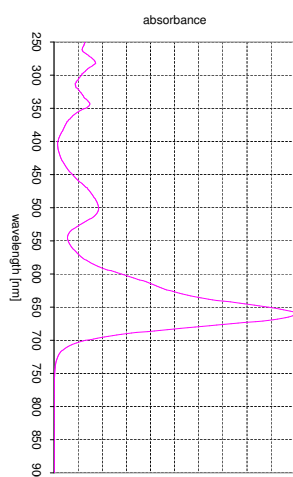
**pH:** 7.5 ± 0.5

**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 EDA-ADP-DYQ-660



absorption spectrum of DYQ 660