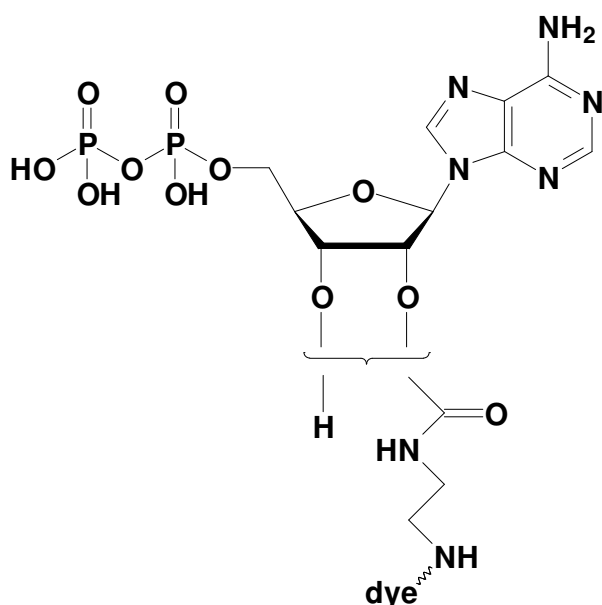


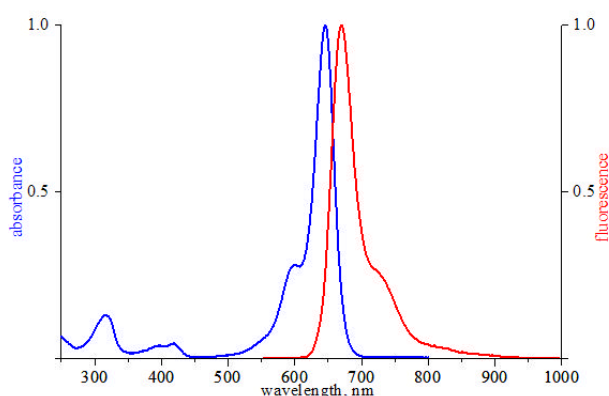
## EDA-ADP - ATTO 647N

2'/3'-O-(2-Aminoethyl-carbamoyl)-Adenosine-5'-diphosphate, labeled with ATTO 647N,  
Triethylammonium salt

Cat. No.	Amount
NU-802-647N	10 $\mu$ l / 1 mM



structural formula of EDA-ADP



excitation and emission spectrum of ATTO 647N

**Cat. No.:** NU-802-647N

**Molecular Formula:** C<sub>13</sub>H<sub>21</sub>N<sub>7</sub>O<sub>11</sub>P<sub>2</sub> - ATTO 647N  
(free acid)

**Molecular Weight:** 1140.29 (free acid)

**Purity:** > 95%, clear aqueous solution, pH 7.5

**Spectroscopic properties:**

$\lambda_{exc}$  644 nm;  $\lambda_{em}$  669 nm;  $\epsilon$  150,000 cm<sup>-1</sup> M<sup>-1</sup>

**Storage conditions:**

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at < -20°C. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

**For research use only!**

**Selected References:**

Hunke *et al.* (2010) The effect of NBD-Cl in nucleotide-binding of the major subunit alpha and B of the motor proteins F1FO ATP synthase and A1AO ATP synthase. *J. Bioenerg. Biomembr.* **41** (1):1.

Luo *et al.* (2008) Crystal structure of the NS3 protease-helicase from dengue virus. *J. Virol.* **82** (1):173.