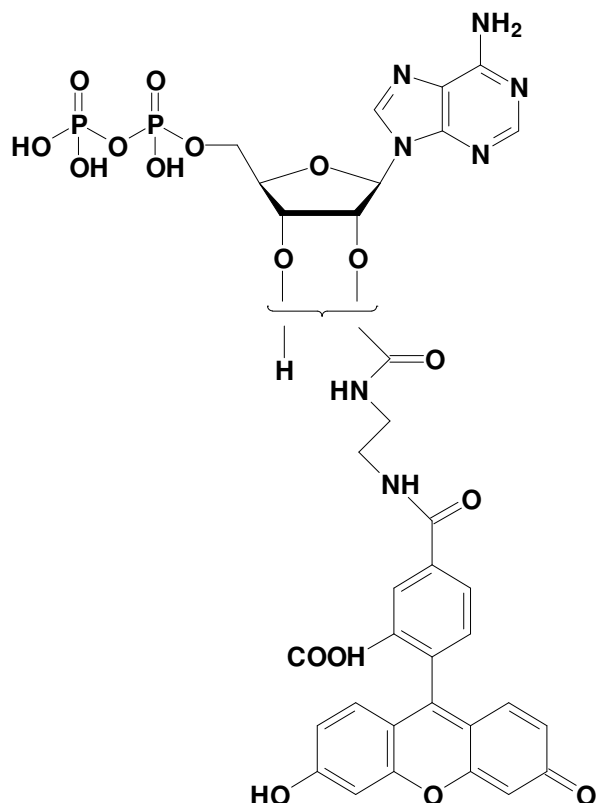


## EDA-ADP - 5 FAM

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

Cat. No.	Amount
NU-802-5FM	160 $\mu$ l (1 mM)



structural formula of EDA-ADP - 5 FAM

**Cat. No.:** NU-802-5FM

**Molecular Formula:** C<sub>34</sub>H<sub>31</sub>N<sub>7</sub>O<sub>17</sub>P<sub>2</sub> (free acid)

**Molecular Weight:** 871.60 (free acid)

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

**Spectroscopic properties:**

$\lambda_{exc}$  492 nm;  $\lambda_{em}$  517 nm;  $\epsilon$  83,000 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!**