



### $\gamma$ -(6-Aminoethyl)-ATP-ATTO-425

$\gamma$ -(6-Aminoethyl)-adenosine-5'-triphosphate, labeled with ATTO 425, Triethylammonium salt

Cat. No.	Amount
NU-833-425	120 $\mu$ l (1 mM)

**For research use only!**

**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>38</sub>H<sub>54</sub>N<sub>7</sub>O<sub>18</sub>P<sub>3</sub> (free acid)

**Molecular Weight:** 989.80 g/mol (free acid)

**Exact Mass:** 989.27 g/mol (free acid)

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

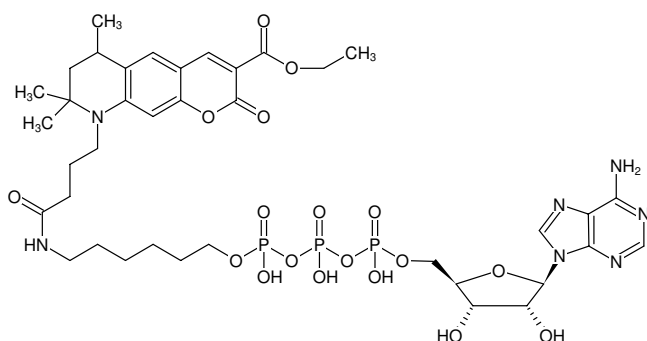
**Form:** solution in water

**Color:** green-yellow

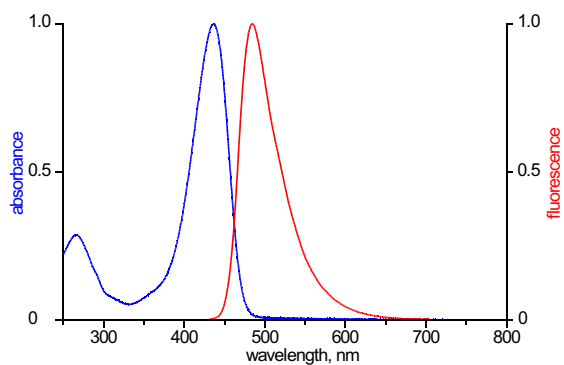
**Concentration:** 1.0 mM - 1.1 mM

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

**Spectroscopic Properties:**  $\lambda_{exc}$  439 nm,  $\lambda_{em}$  485 nm,  $\epsilon$  45.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)



Structural formula of  $\gamma$ -(6-Aminoethyl)-ATP-ATTO-425



excitation and emission spectrum of ATTO 425