

**EDA-ATP - ATTO-540Q**

2'/3'-O-(2-Aminoethyl-carbamoyl)-Adenosine-5'-triphosphate,  
labeled with ATTO 540Q, Triethylammonium salt

Cat. No.	Amount
NU-808-540Q	40 µl (1 mM)

**For research use only!**

**Shipping:** shipped on blue ice

**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>13</sub>H<sub>22</sub>N<sub>7</sub>O<sub>14</sub>P<sub>3</sub> - ATTO 540Q (free acid)

**Molecular Weight:** 1133.27 g/mol (free acid)

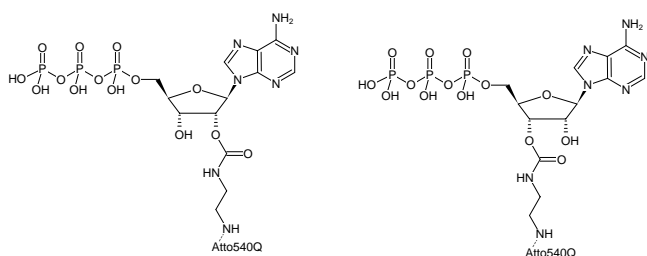
**Purity:** ≥ 95 % (HPLC)

**Form:** clear aqueous solution

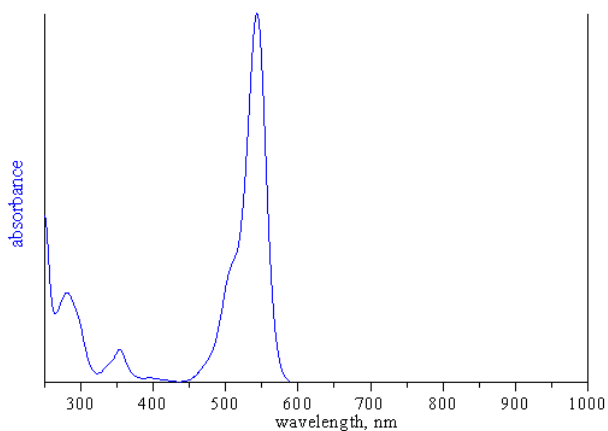
**Concentration:** 1 mM - 1.1 mM

**pH:** 7.5 ± 0.5

**Spectroscopic Properties:** λ<sub>exc</sub> 543 nm, ε 105.0 L mmol<sup>-1</sup> cm<sup>-1</sup>  
(Tris-HCl pH 7.5)



Structural formula of EDA-ATP - ATTO-540Q



absorption spectrum of ATTO 540Q