

**EDA-ATPyS-ATTO-740**

2'/3'-O-(2-Aminoethyl-carbamoyl)-adenosine-5'-(γ-thio)-triphosphate, labeled with ATTO 740, Triethylammonium salt

Cat. No.	Amount
NU-1609-740	50 µl (1 mM)

**For research use only!**

**Shipping:** shipped on dry ice

**Storage Conditions:** store at -20 °C

**Additional Storage Conditions:** long time storage at pH 7.5

**Shelf Life:** 6 months after date of delivery

**Molecular Formula:** C<sub>13</sub>H<sub>22</sub>N<sub>7</sub>O<sub>13</sub>P<sub>3</sub>S - ATTO 740 (free acid)

**Molecular Weight:** 1058.33 g/mol (free acid)

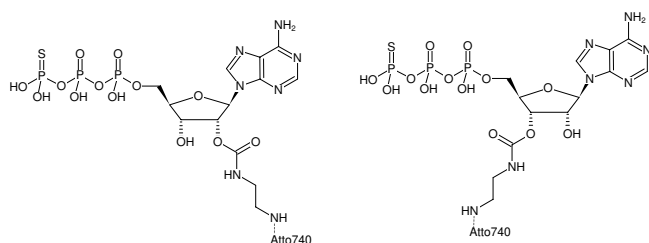
**Purity:** ≥ 90 % (HPLC), contains approx. 5 % EDA-ADP-dye

**Form:** solution in water

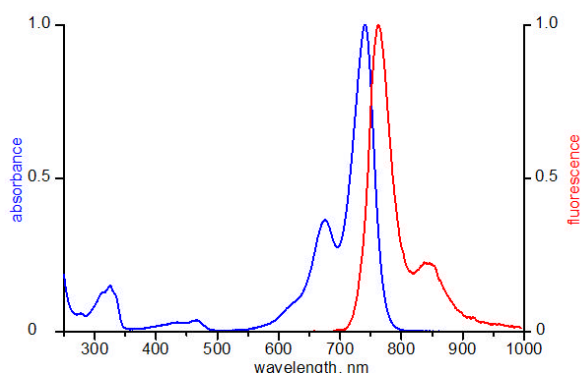
**Concentration:** 1.0 mM - 1.1 mM

**pH:** 7.5 ±0.5

**Spectroscopic Properties:** λ<sub>exc</sub> 743 nm, λ<sub>em</sub> 763 nm, ε 120.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)



Structural formula of EDA-ATPyS-ATTO-740



excitation and emission spectrum of ATTO 740