



## EDA-ATPyS-6-FAM

2'/3'-O-(2-Aminoethyl-carbamoyl)-adenosine-5'-(γ-thio)-triphosphate, labeled with 6 FAM, Triethylammonium salt

Cat. No.	Amount
NU-1609-6FM	50 µl (1 mM)

**For general laboratory use.**

**Shipping:** shipped on dry ice

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

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

**Molecular Formula:** C<sub>34</sub>H<sub>32</sub>N<sub>7</sub>O<sub>19</sub>P<sub>3</sub>S (free acid)

**Molecular Weight:** 967.64 g/mol (free acid)

**Purity:** ≥ 90 % (HPLC)

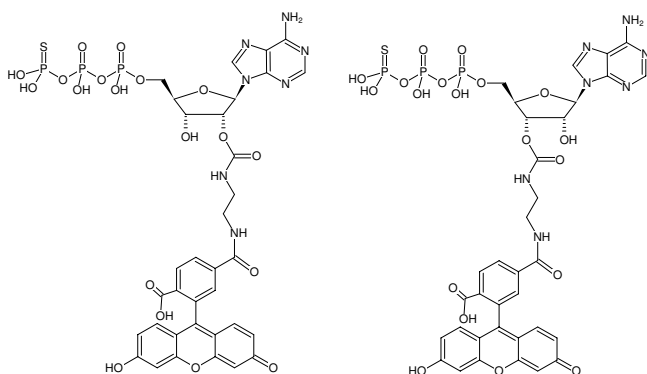
**Form:** solution in water

**Color:** yellow

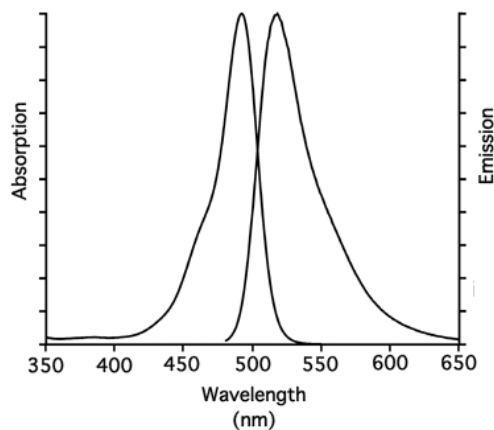
**Concentration:** 1.0 mM - 1.1 mM

**pH:** 7.5 ±0.5

**Spectroscopic Properties:** λ<sub>exc</sub> 492 nm, λ<sub>em</sub> 517 nm, ε 83.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)



Structural formula of EDA-ATPyS-6-FAM



excitation and emission spectrum of 6-FAM