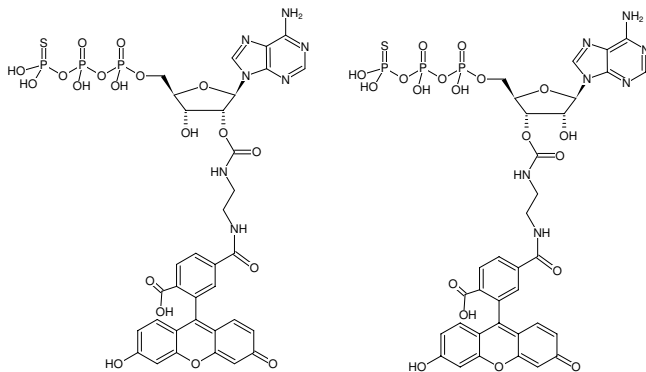




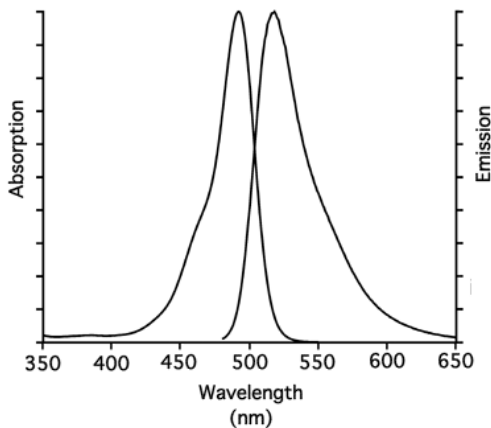
EDA-ATPyS-6-FAM

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)



Structural formula of EDA-ATPyS-6-FAM



excitation and emission spectrum of 6-FAM

For research use only!

Shipping: shipped on dry ice

Storage Conditions: store at -20 °C

Shelf Life: 6 months after date of delivery

Molecular Formula: C₃₄H₃₂N₇O₁₉P₃S (free acid)

Molecular Weight: 967.64 g/mol (free acid)

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

Form: solution in water

Color: yellow

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 492 nm, λ_{em} 517 nm, ε 83.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)

Please note: For reasons of stability, please make sure that the pH value of a solution of this product never drops below 7.0.

The limited stability of labeled ATPγS and GTPγS nucleotides makes it impossible to keep stocks of them and to offer them as catalog products that are available immediately. Therefore, any labeled ATPγS or GTPγS is freshly prepared per order. For this reason the price is higher than for stable nucleotide-dye conjugates which can be synthesized and kept in stock.