



## EDA-ATPyS-Cy5

2'/3'-O-(2-Aminoethyl-carbamoyl)-adenosine-5'-(γ-thio)-triphosphate, labeled with Cy5, Triethylammonium salt

| Cat. No.    | Amount       |
|-------------|--------------|
| NU-1609-CY5 | 50 µl (1 mM) |

**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:** C<sub>46</sub>H<sub>60</sub>N<sub>9</sub>O<sub>20</sub>P<sub>3</sub>S<sub>3</sub> (free acid)

**Molecular Weight:** 1248.13 g/mol (free acid)

**Exact Mass:** 1247.23 g/mol (free acid)

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

**Form:** solution in water

**Color:** blue

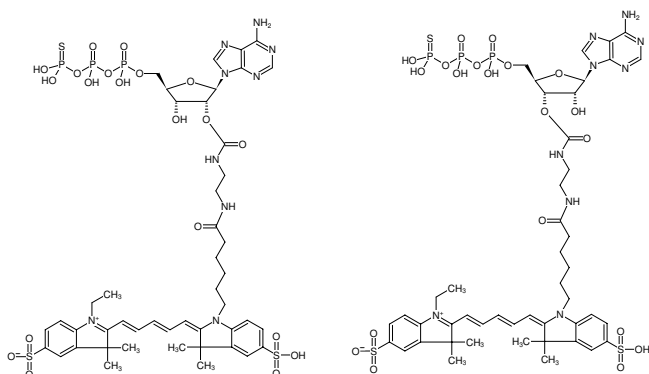
**Concentration:** 1.0 mM - 1.1 mM

**pH:** 7.5 ± 0.5

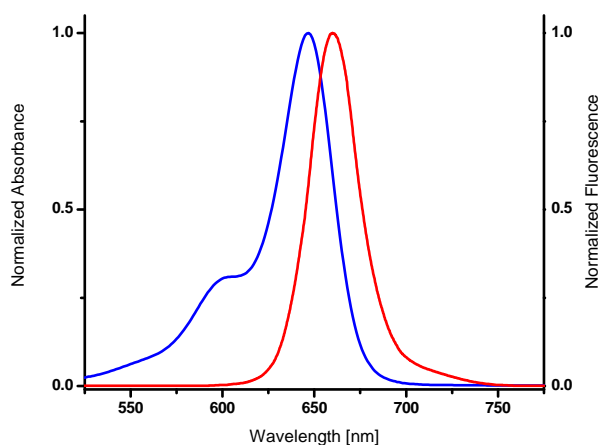
**Spectroscopic Properties:** λ<sub>exc</sub> 649 nm, λ<sub>em</sub> 670 nm, ε 250.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)

**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 ATPyS and GTPyS nucleotides makes it impossible to keep stocks of them and to offer them as catalog products that are available immediately. Therefore, any labeled ATPyS or GTPyS 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.



Structural formula of EDA-ATPyS-Cy5



excitation and emission spectrum of Cy5