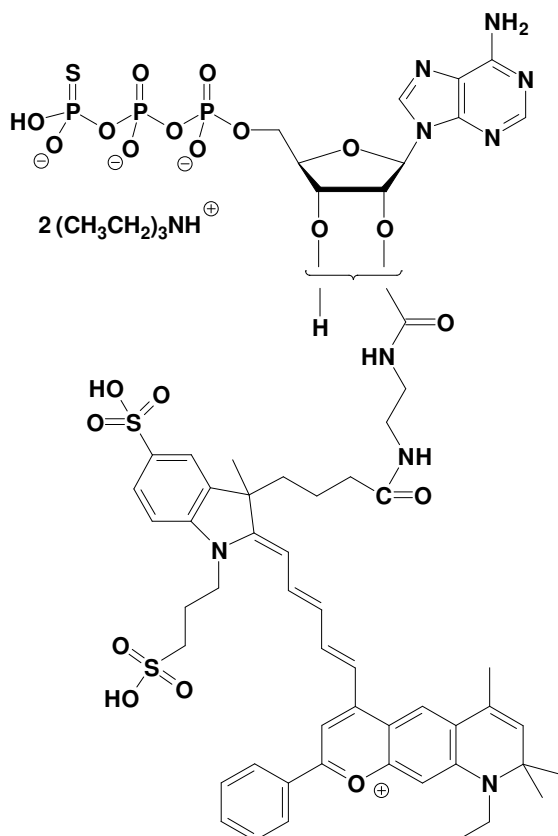


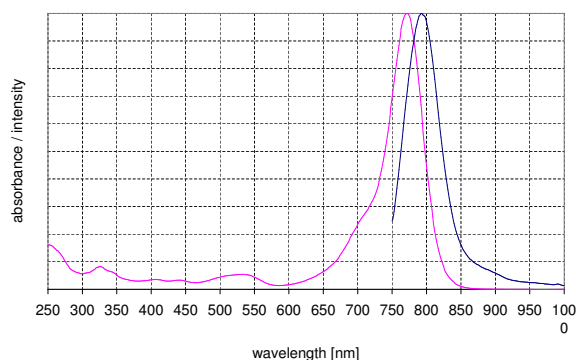
## EDA-ATP $\gamma$ S - DY 776

2'/3'-O-(2-Aminoethyl-carbamoyl)-Adenosine-5'-( $\gamma$ -thio)-triphosphate, labeled with DY 776, Triethylammonium salt

Cat. No.	Amount
NU-1609-776	50 $\mu$ l / 1 mM



structural formula of EDA-ATP $\gamma$ S - DY 776



excitation and emission spectrum of DY 776

**Cat. No.:** NU-1609-776

**Molecular Formula:** C<sub>57</sub>H<sub>66</sub>N<sub>9</sub>O<sub>21</sub>P<sub>3</sub>S<sub>3</sub> (Anion)

**Molecular Weight:** 1402.31 (Anion)

**Purity:** > 95%, clear aqueous solution, pH 7.5

**Spectroscopic properties:**

$\lambda_{exc}$  771 nm;  $\lambda_{em}$  801 nm;  $\epsilon$  120,000 cm<sup>-1</sup> M<sup>-1</sup>

**Storage conditions:**

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at < -20°C. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

**For research use only!**