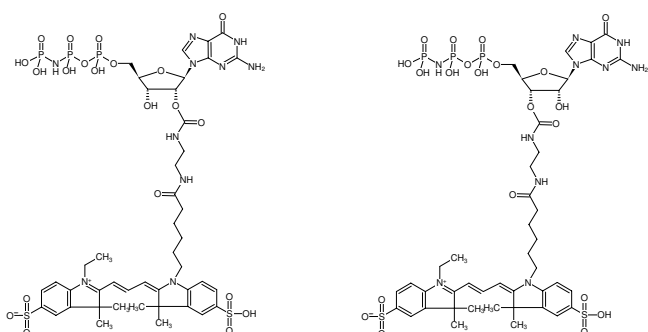




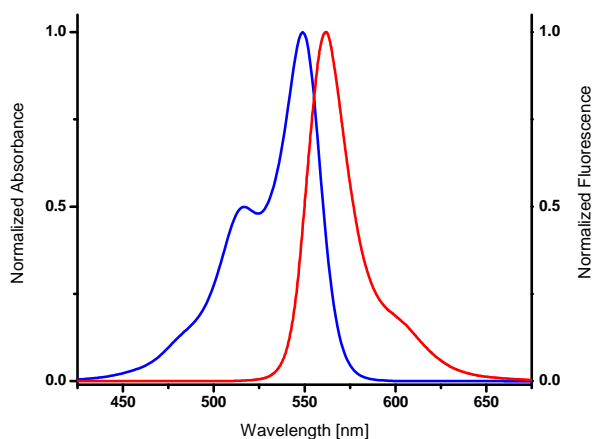
## EDA-GppNHp (EDA-GMPPNP)-Cy3

2'/3'-O-(2-Aminoethyl-carbamoyl)-Guanosine-5'-[( $\beta,\gamma$ -imido)] triphosphate, labeled with Cy3, Triethylammonium salt

Cat. No.	Amount
NU-860-CY3	50 $\mu$ l (1 mM)



Structural formula of EDA-GppNHp (EDA-GMPPNP)-Cy3



excitation and emission spectrum of Cy3

**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>44</sub>H<sub>59</sub>N<sub>10</sub>O<sub>21</sub>P<sub>3</sub>S<sub>2</sub> (free acid)

**Molecular Weight:** 1221.05 g/mol (free acid)

**Exact Mass:** 1220.25 g/mol (free acid)

**Purity:**  $\geq$  95 % (HPLC)

**Form:** solution in water

**Color:** pink

**Concentration:** 1.0 mM - 1.1 mM

**pH:** 7.5  $\pm$  0.5

**Spectroscopic Properties:**  $\lambda_{exc}$  550 nm,  $\lambda_{em}$  570 nm,  $\epsilon$  150.0 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. This can be achieved by dissolving the nucleotide in a buffer of your choice (50 - 100 mM, pH 7 - 10). Dissolve and adjust concentration photometrically.