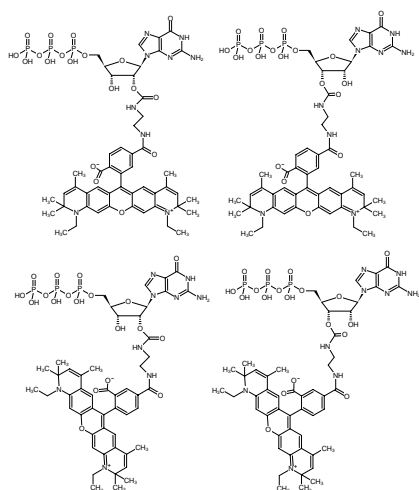




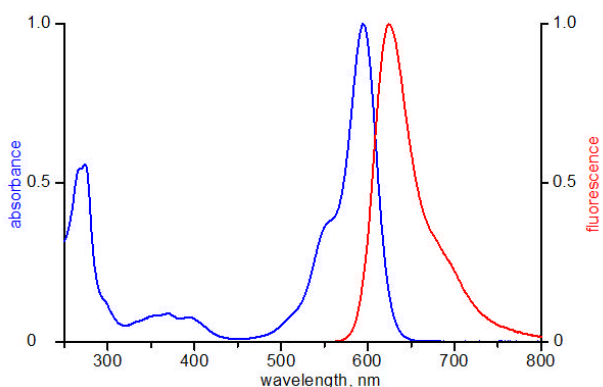
EDA-GTP - ATTO-590

2'/3'-O-(2-Aminoethyl-carbamoyl)-Guanosine-5'-triphosphate, labeled with ATTO 590, Triethylammonium salt

Cat. No.	Amount
NU-820-590	120 µl (1 mM)



Structural formula of EDA-GTP - ATTO-590



excitation and emission spectrum of ATTO 590

For research use only!

Shipping: shipped on blue ice

Storage Conditions: store at -20 °C

Short term exposure (up to 1 week cumulative) to ambient temperature possible.

Shelf Life: 12 months after date of delivery

Molecular Formula: C₅₀H₅₈N₉O₁₉P₃ (free acid)

Molecular Weight: 1181.98 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: clear aqueous solution

Concentration: 1.0 mM - 11 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 593 nm, λ_{em} 622 nm, ε 120.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)

Applications:

Myosin labelling and fluorometric binding studies^[1]

Nucleotide exchange at myosin^[2]

Selected References:

- [1] Oiwa *et al.* (2003) Myosin regulatory light chain phosphorylation and strain modulate adenosin diphosphate release from smooth muscle myosin. *Bioph. J.* **84**:634.
- [2] Shirakava *et al.* (2000) Measurement of nucleotide exchange rate constants in single rabbit soleus myofibrils during shortening and lengthening using a fluorescent ATP. *Biophys. J.* **78**:218.