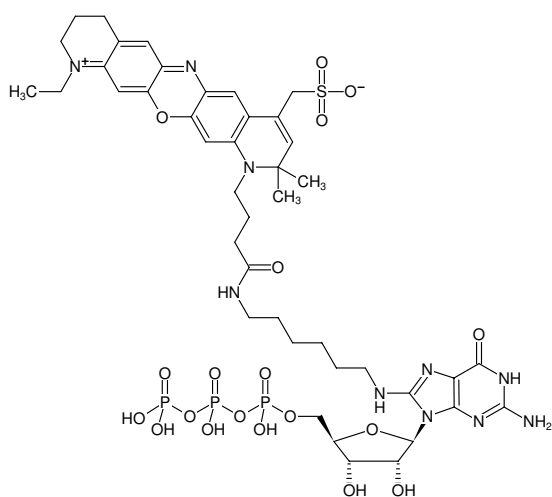




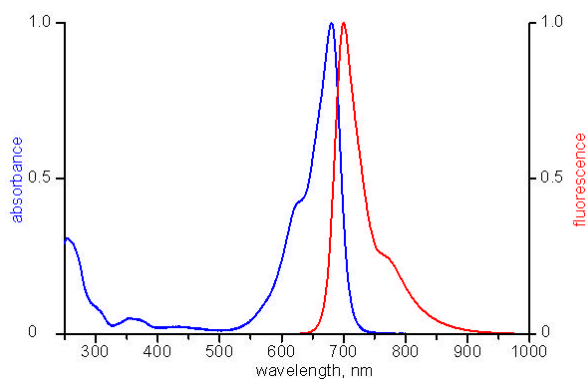
8-(6-Aminoethyl)-amino-GTP-ATTO-680

8-(6-Aminoethyl)-amino-guanosine-5'-triphosphate, labeled with ATTO 680, Triethylammonium salt

Cat. No.	Amount
NU-830-680	80 µl (1 mM)



Structural formula of 8-(6-Aminoethyl)-amino-GTP-ATTO-680



excitation and emission spectrum of ATTO 680

For general laboratory use.

Shipping: shipped on gel packs

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: C₄₃H₅₉N₁₀O₁₉P₃S (free acid)

Molecular Weight: 1144.98 g/mol (free acid)

Exact Mass: 1144.29 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: solution in water

Color: blue

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 681 nm, λ_{em} 698 nm, ε 125.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)

Selected References:

Feibelman *et al.* (2018) Identification of small molecule inhibitors of the Chikungunya virus nsP1 RNA capping enzyme. *Antiviral Res.* **154**:124.

Bullard-Feibelman *et al.* (2016) A Sensitive and Robust High-Throughput Screening Assay for Inhibitors of the Chikungunya Virus nsP1 Capping Enzyme. *PLoS One* **11** (7):e0158923.

Geiss *et al.* (2011) A High-Throughput Screening Assay for the Identification of Flavivirus NS5 Capping Enzyme GTP-Binding Inhibitors: Implications for Antiviral Drug Development. *Journal of Biomolecular Screening* **16**(8):852.