

**EDA-GTP-ATTO-550**

2'/3'-O-(2-Aminoethyl-carbamoyl)-Guanosine-5'-triphosphate, labeled with ATTO 550, Triethylammonium salt

Cat. No.	Amount
NU-820-550	160 µl (0,5 mM)

**For research use only!**

**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:** C<sub>50</sub>H<sub>63</sub>N<sub>10</sub>O<sub>18</sub>P<sub>3</sub> (free acid)

**Molecular Weight:** 1185.03 g/mol (free acid)

**Exact Mass:** 1184.35 g/mol (free acid)

**Purity:** ≥ 95 % (HPLC)

**Form:** solution in water, containing 10% DMSO

**Color:** red-violet

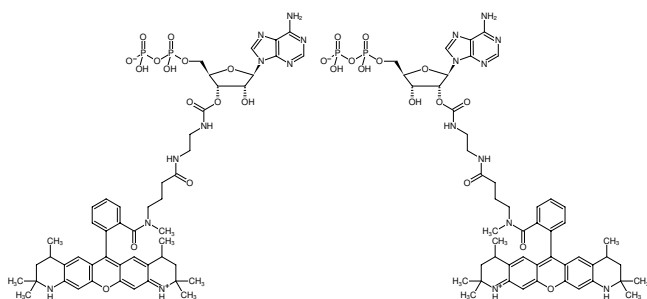
**Concentration:** 0.50 mM - 0.55 mM

**pH:** 7.5 ± 0.5

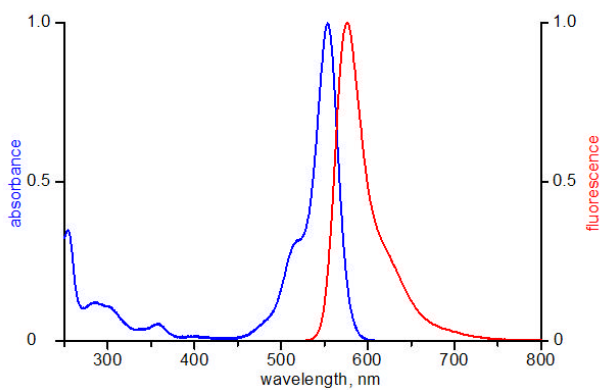
**Spectroscopic Properties:** λ<sub>exc</sub> 554 nm, λ<sub>em</sub> 576 nm, ε 120.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)

**Selected References:**

Oreb *et al.* (2011) Substrate binding disrupts dimerization and induces nucleotide exchange of the chloroplast GTPase Toc33. *J. Biol. Chem.* **436** (2):31.



Structural formula of EDA-GTP-ATTO-550



excitation and emission spectrum of ATTO 550