



## EDA-GTPyS-ATTO-680

2'/3'-O-(2-Aminoethyl-carbamoyl)-guanosine-5'-(γ-thio)-triphosphate, labeled with ATTO 680, Triethylammonium salt

Cat. No.	Amount
NU-1610-680	50 µl (1 mM)

**For research use only!**

**Shipping:** shipped on dry ice

**Storage Conditions:** store at -20 °C

**Shelf Life:** 6 months after date of delivery

**Molecular Formula:** C<sub>40</sub>H<sub>51</sub>N<sub>10</sub>O<sub>19</sub>P<sub>3</sub>S<sub>2</sub> (free acid)

**Molecular Weight:** 1132.94 g/mol (free acid)

**Exact Mass:** 1132.20 g/mol (free acid)

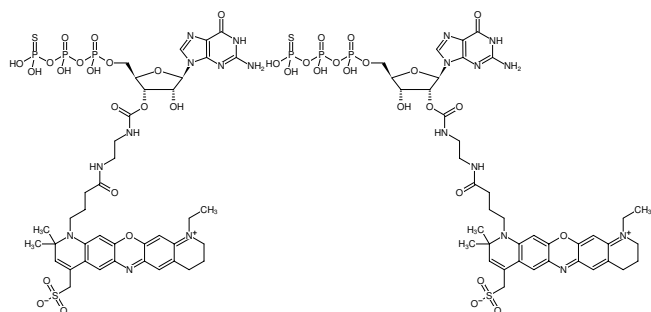
**Purity:** ≥ 90 % (HPLC), contains approx. 5 % EDA-GDP-dye

**Form:** solution in water

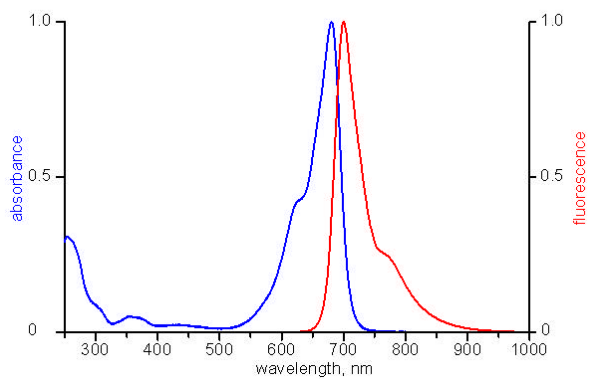
**Concentration:** 1.0 mM - 1.1 mM

**pH:** 7.5 ± 0.5

**Spectroscopic Properties:** λ<sub>exc</sub> 681 nm, λ<sub>em</sub> 698 nm, ε 125.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)



Structural formula of EDA-GTPyS-ATTO-680



excitation and emission spectrum of ATTO 680