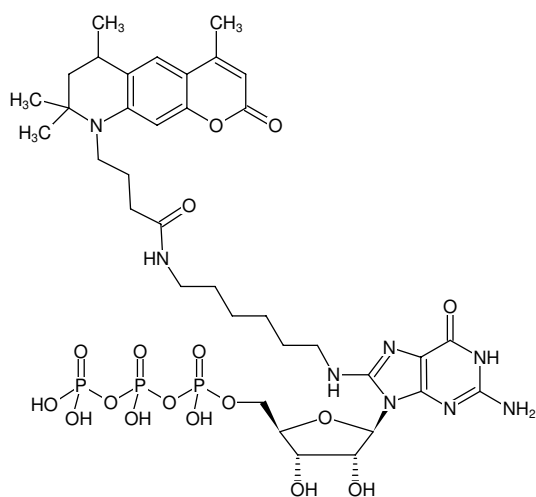




## 8-(6-Aminoethyl)-amino-GTP-ATTO-390

8-(6-Aminoethyl)-amino-guanosine-5'-triphosphate, labeled with ATTO 390 Triethylammonium salt

Cat. No.	Amount
NU-830-390	120 µl (1 mM)



Structural formula of 8-(6-Aminoethyl)-amino-GTP-ATTO-390

**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:** C<sub>36</sub>H<sub>53</sub>N<sub>8</sub>O<sub>17</sub>P<sub>3</sub> (free acid)

**Molecular Weight:** 962.78 g/mol (free acid)

**Exact Mass:** 962.27 g/mol (free acid)

**Purity:** ≥ 95 % (HPLC)

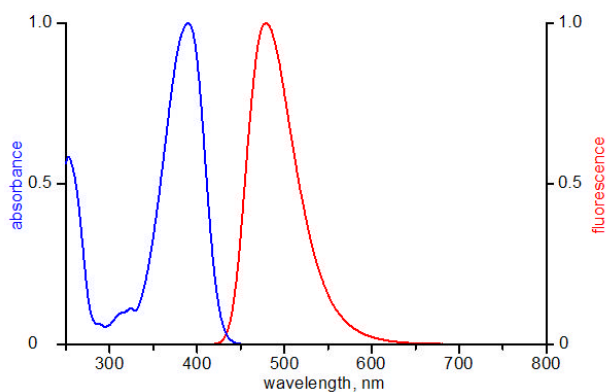
**Form:** solution in water

**Color:** slightly yellow

**Concentration:** 1.0 mM - 1.1 mM

**pH:** 7.5 ± 0.5

**Spectroscopic Properties:** λ<sub>exc</sub> 390 nm, λ<sub>em</sub> 476 nm, ε 24.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)



excitation and emission spectrum of ATTO 390