



## 8-(6-Aminoethyl)-amino-GMP-ATTO-590

8-(6-Aminoethyl)-amino-guanosine-5'-monophosphate, labeled with ATTO 590

Cat. No.	Amount
NU-829-590	120 µl (1 mM)

**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:** C<sub>53</sub>H<sub>64</sub>N<sub>9</sub>O<sub>12</sub>P (free acid)

**Molecular Weight:** 1050.11 g/mol (free acid)

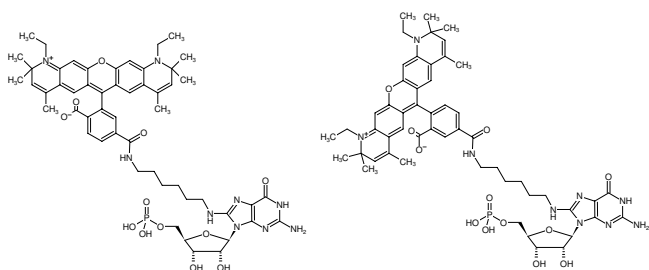
**Purity:** ≥ 95 % (HPLC)

**Form:** solution in water

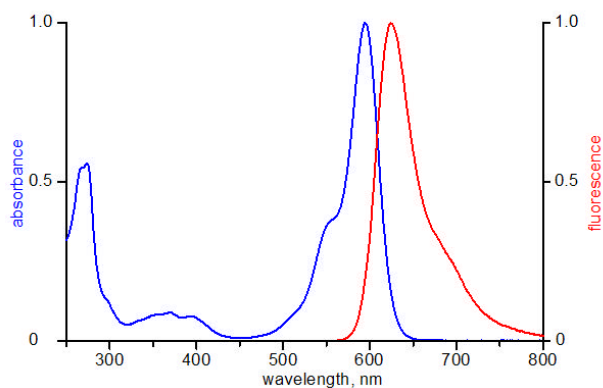
**Concentration:** 1.0 mM - 1.1 mM

**pH:** 7.5 ± 0.5

**Spectroscopic Properties:** λ<sub>exc</sub> 593 nm, λ<sub>em</sub> 622 nm, ε 120.0 L mol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)



Structural formula of 8-(6-Aminoethyl)-amino-GMP-ATTO-590



excitation and emission spectrum of ATTO 590