



### 8-(6-Aminoethyl)-amino-GMP-ATTO-612Q

8-(6-Aminoethyl)-amino-guanosine-5'-monophosphate, labeled with ATTO 612Q, Triethylammonium salt

Cat. No.	Amount
NU-829-612Q	80 µl (1 mM)

For general laboratory use.

Please centrifuge briefly before opening (volume ≤2 ml).

**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>16</sub>H<sub>28</sub>N<sub>7</sub>O<sub>8</sub>P - ATTO 612Q (free acid)

**Molecular Weight:** 1149.41 g/mol (free acid)

**Purity:** ≥ 95 % (HPLC)

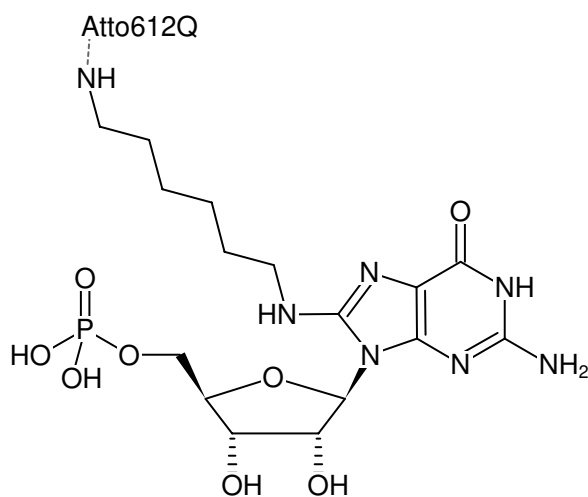
**Form:** solution in water

**Color:** blue

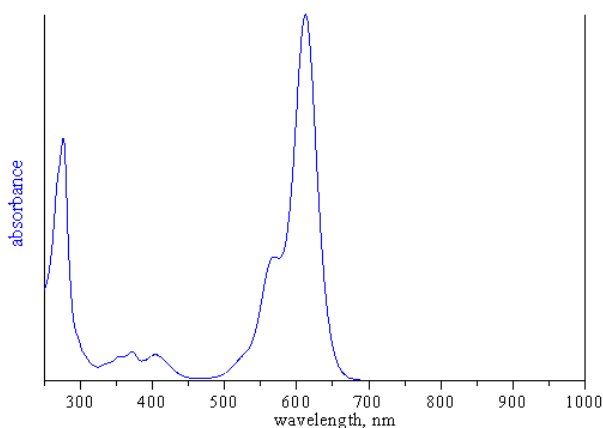
**Concentration:** 1.0 mM - 1.1 mM

**pH:** 7.5 ±0.5

**Spectroscopic Properties:** λ<sub>exc</sub> 615 nm, ε 115.0 L mol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)



Structural formula of 8-(6-Aminoethyl)-amino-GMP-ATTO-612Q



absorption spectrum of ATTO 612Q