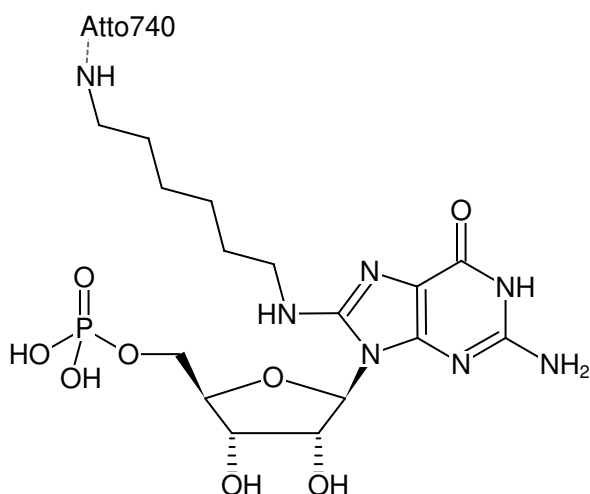




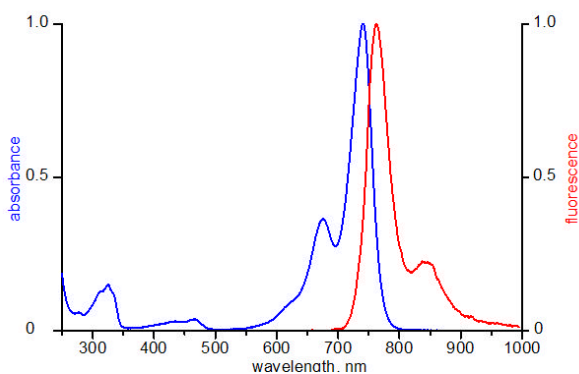
## 8-(6-Aminoethyl)-amino-GMP-ATTO-740

8-(6-Aminoethyl)-amino-guanosine-5'-monophosphate, labeled with ATTO 740, Triethylammonium salt

Cat. No.	Amount
NU-829-740	80 µl (1 mM)



Structural formula of 8-(6-Aminoethyl)-amino-GMP-ATTO-740



excitation and emission spectrum of ATTO 740

**For general laboratory use.**

**Shipping:** shipped on gel packs

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

**Additional Storage Conditions:** long time storage at pH 7.5

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 740 (free acid)

**Molecular Weight:** 926.41 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> 743 nm, λ<sub>em</sub> 763 nm, ε 120.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)

**Please note:** The dye decomposes at pH-levels greater 8.0. The recommended pH for a solution of this product is 7.5.