



### N<sup>6</sup>-(6-Aminoethyl)-ATP-ATTO-643

N<sup>6</sup>-(6-Aminoethyl)-adenosine-5'-triphosphate, labeled with ATTO 643, Triethylammonium salt

Cat. No.	Amount
NU-805-643	120 µl (1 mM)

**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>16</sub>H<sub>28</sub>N<sub>6</sub>O<sub>13</sub>P<sub>3</sub> - ATTO 643 (free acid)

**Molecular Weight:** 1423.45 g/mol - ATTO 643 (free acid)

**Purity:** ≥ 95 % (HPLC)

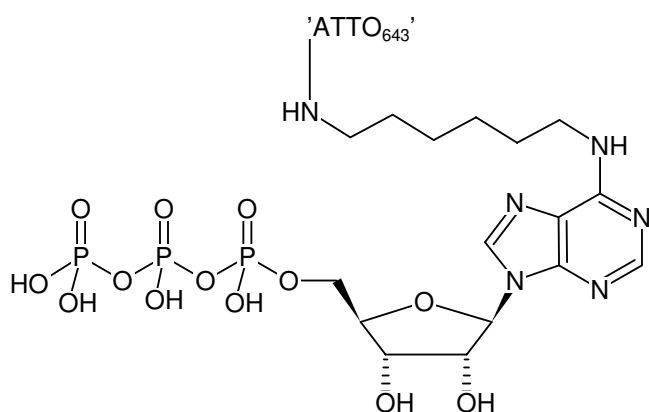
**Form:** solution in water

**Color:** blue

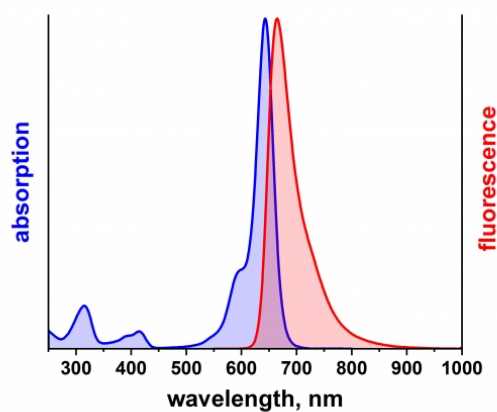
**Concentration:** 1.0 mM - 1.1 mM

**pH:** 7.5 ± 0.5

**Spectroscopic Properties:** λ<sub>exc</sub> 643 nm, λ<sub>em</sub> 665 nm, ε 150.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)



Structural formula of N<sup>6</sup>-(6-Aminoethyl)-ATP-ATTO-643



excitation and emission spectrum of ATTO 643