



### N<sup>6</sup>-(6-Aminoheptyl)-dATP-ATTO-580Q

N<sup>6</sup>-(6-Aminoheptyl)-2'-deoxyadenosine-5'-triphosphate, labeled with ATTO 580Q, Triethylammonium salt

Cat. No.	Amount
NU-835-580Q	40 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>29</sub>N<sub>6</sub>O<sub>12</sub>P<sub>3</sub> - ATTO 580Q (free acid)

**Molecular Weight:** 1266.36 g/mol (free acid)

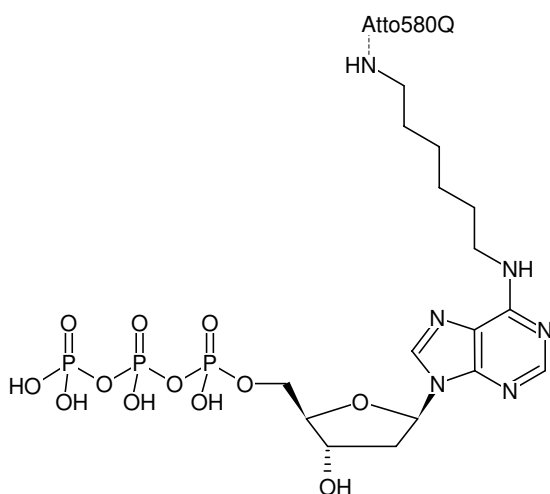
**Purity:** ≥ 95 % (HPLC)

**Form:** solution in water

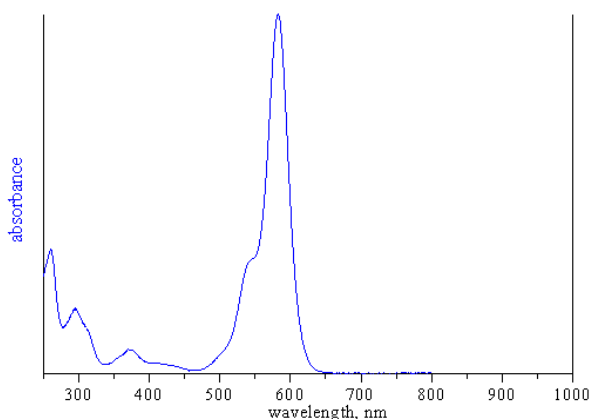
**Concentration:** 1.0 mM - 1.1 mM

**pH:** 7.5 ± 0.5

**Spectroscopic Properties:** λ<sub>exc</sub> 587 nm, ε 110.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)



Structural formula of N<sup>6</sup>-(6-Aminoheptyl)-dATP-ATTO-580Q



absorption spectrum of ATTO 580Q