



### 8-[(6-Amino)hexyl]-amino-cAMP - ATTO-540Q

8-[(6-Amino)hexyl]-amino-adenosine-3',5'-cyclic monophosphate, labeled with ATTO 540Q, Triethylammonium salt

Cat. No.	Amount
NU-851-540Q	40 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>16</sub>H<sub>26</sub>N<sub>7</sub>O<sub>6</sub>P - ATTO 540Q (free acid)

**Molecular Weight:** 983.39 g/mol (free acid)

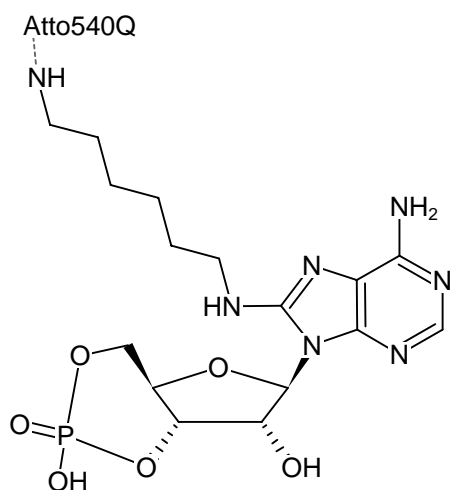
**Purity:** ≥ 95 % (HPLC)

**Form:** clear aqueous solution, pH 7.5 ± 0.5

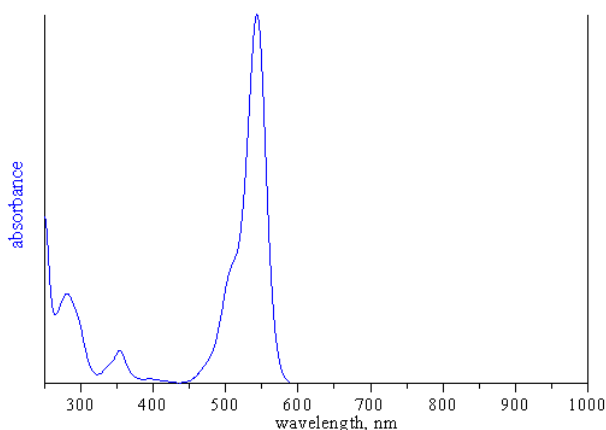
**Concentration:** 1 mM - 1.1 mM

**pH:** 7.5 ± 0.5

**Spectroscopic Properties:** λ<sub>exc</sub> 543 nm, ε 105.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)



Structural formula of 8-[(6-Amino)hexyl]-amino-cAMP - ATTO-540Q



absorption spectrum of ATTO 540Q