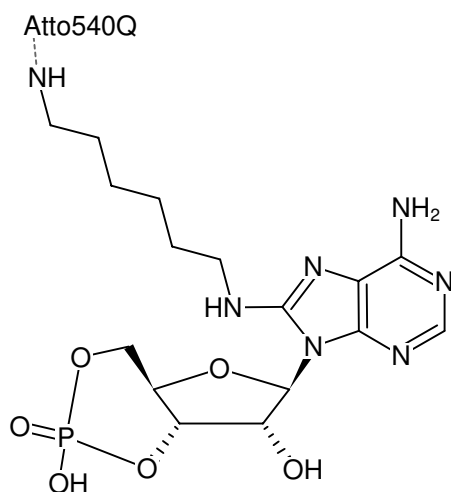




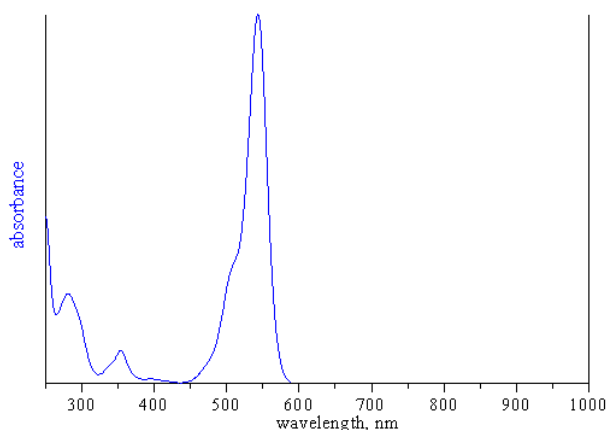
8-(6-Aminoheptyl)-amino-cAMP-ATTO-540Q

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

Cat. No.	Amount
NU-851-540Q	40 µl (1 mM)



Structural formula of 8-(6-Aminoheptyl)-amino-cAMP-ATTO-540Q



absorption spectrum of ATTO 540Q

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: C₁₆H₂₆N₇O₆P - ATTO 540Q (free acid)

Molecular Weight: 983.39 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: clear aqueous solution

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 543 nm, ε 105.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)