


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. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

Shelf Life: 12 months

Molecular Formula: C₁₆H₂₆N₇O₆P - ATTO 540Q (free acid)

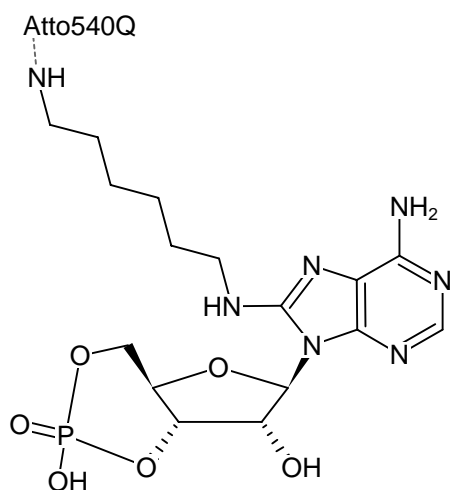
Molecular Weight: 983.39 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

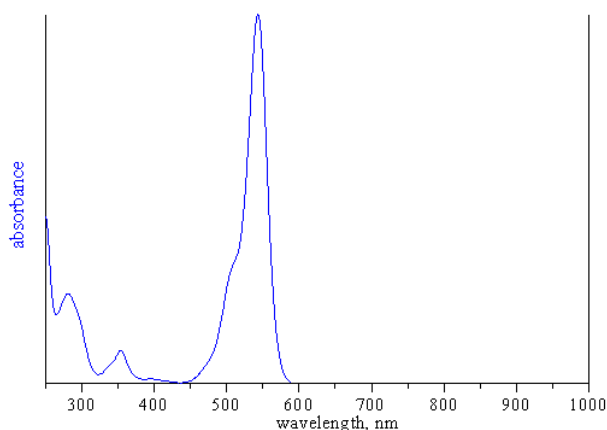
Form: clear aqueous solution, pH 7.5 ± 0.5

Concentration: 1 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 542 nm; ε 105.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)


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



absorption spectrum of ATTO 540Q