



N⁶-(6-Aminohexyl)-ADP-ATTO-643

N⁶-(6-Aminohexyl)-adenosine-5'-diphosphate, labeled with ATTO 643, Triethylammonium salt

Cat. No.	Amount
NU-247-643	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: C₁₆H₂₇N₆O₁₀P₂ - ATTO 643 (free acid)

Molecular Weight: 1344.7 g/mol

Purity: ≥ 95 % (HPLC)

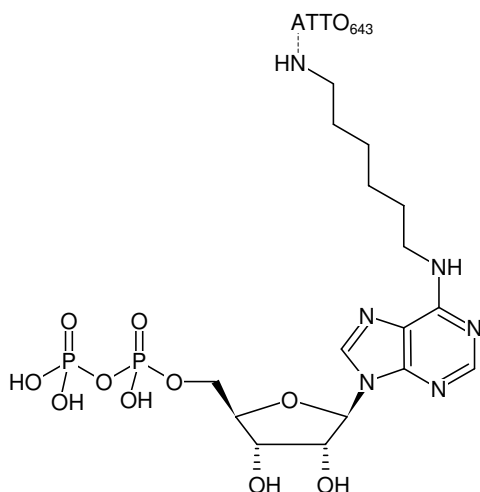
Form: solution in water

Color: blue

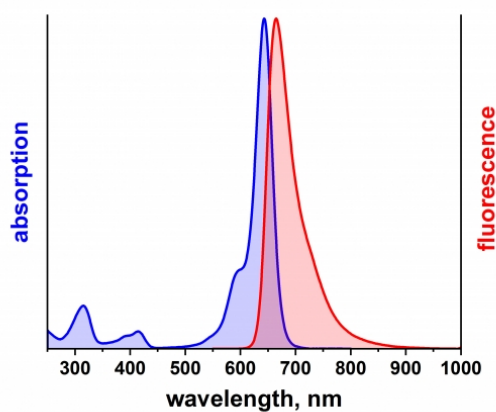
Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 643 nm, λ_{em} 665 nm, ε 150.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



Structural formula of N⁶-(6-Aminohexyl)-ADP-ATTO-643



excitation and emission spectrum of ATTO 643