



N⁶-(6-Aminohexyl)-ATP-ATTO-700

N⁶-(6-Aminohexyl)-adenosine-5'-triphosphate, labeled with ATTO 700, Triethylammonium salt

| Cat. No. | Amount |
|------------|--------------|
| NU-805-700 | 80 µl (1 mM) |

For general laboratory use.

Please centrifuge briefly before opening (volume ≤2 ml).

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: C₄₆H₆₂N₉O₁₈P₃S (free acid)

Molecular Weight: 1154.03 g/mol (free acid)

Exact Mass: 1153.31 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

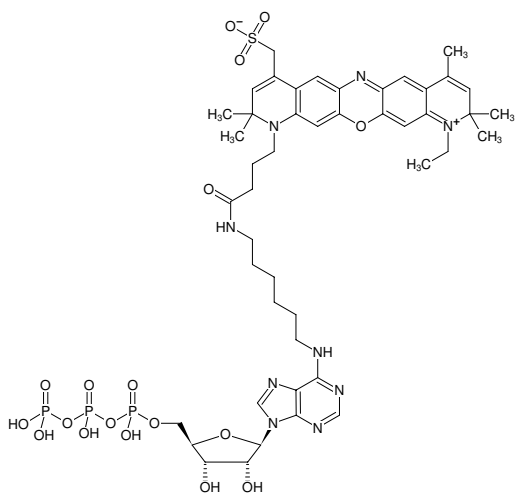
Form: solution in water

Color: blue

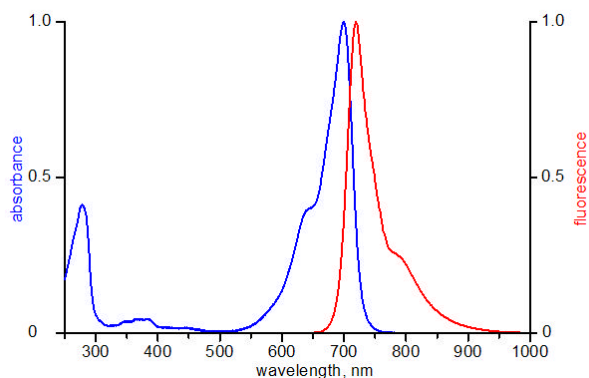
Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ±0.5

Spectroscopic Properties: λ_{exc} 700 nm, λ_{em} 716 nm, ε 120.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



Structural formula of N⁶-(6-Aminohexyl)-ATP-ATTO-700



excitation and emission spectrum of ATTO 700