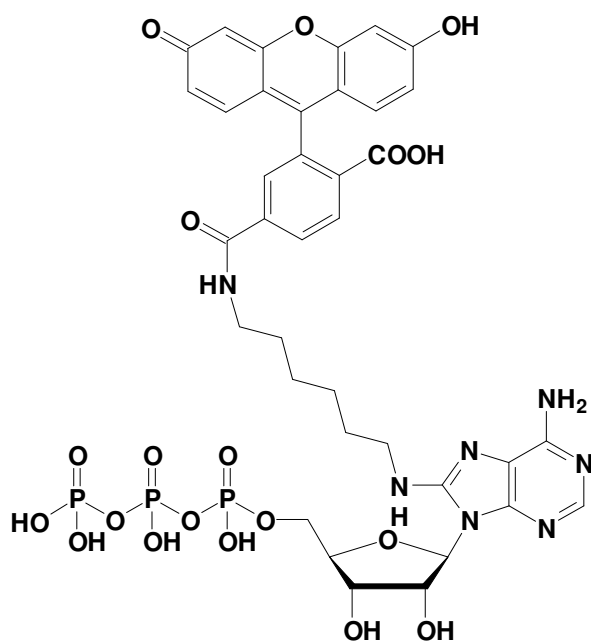


8-[(6-Amino)hexyl]-amino-ATP - 6 FAM

8-[(6-Amino)hexyl]-amino-adenosine-5'-triphosphate, labeled with 6 FAM,
 Triethylammonium salt

Cat. No.	Amount
NU-807-6FM	160 μ l (1 mM)



structural formula of 8-[(6-Amino)hexyl]-amino-ATP - 6 FAM

Cat. No.: NU-807-6FM

Molecular Formula: C₃₇H₄₀N₇O₁₉P₃ (free acid)

Molecular Weight: 979.68 (free acid)

Purity: > 95%, clear aqueous solution, pH 7.5

Spectroscopic properties:

λ_{exc} 492 nm; λ_{em} 517 nm; ϵ 83,000 cm⁻¹ M⁻¹

Storage conditions:

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at < -20°C. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

For research use only!