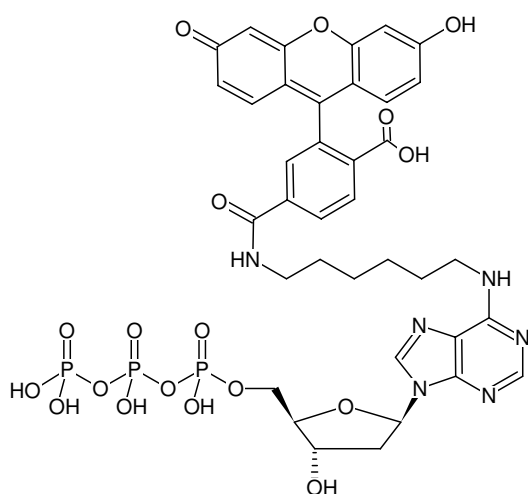




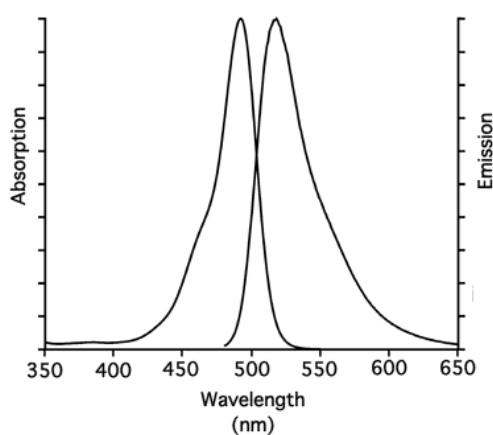
N⁶-(6-Aminoheptyl)-dATP-6-FAM

N⁶-(6-Aminoheptyl)-2'-deoxyadenosine-5'-triphosphate, labeled with 6 FAM, Triethylammonium salt

| Cat. No. | Amount |
|------------|---------------|
| NU-835-6FM | 160 µl (1 mM) |



Structural formula of N⁶-(6-Aminoheptyl)-dATP-6-FAM



excitation and emission spectrum of 6-FAM

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: C₃₇H₃₉N₆O₁₈P₃ (free acid)

Molecular Weight: 948.66 g/mol (free acid)

Exact Mass: 948.15 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: solution in water

Color: yellow

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 492 nm, λ_{em} 517 nm, ε 83.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)

Selected References:

Bianchi *et al.* (2000) PrimPol bypasses UV photoproducts during eukaryotic chromosomal DNA replication. *Mol. Cell* **52** (4):566.