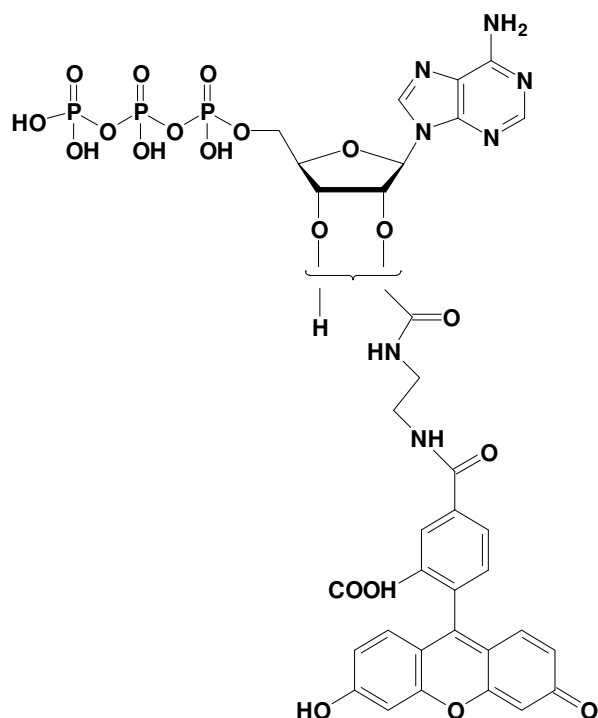


EDA-ATP - 5 FAM

2'/3'-O-(2-Aminoethyl-carbamoyl)-Adenosine-5'-triphosphate, labeled with 5 FAM,
 Triethylammonium salt

| Cat. No. | Amount |
|------------|--------------------|
| NU-808-5FM | 160 μ l (1 mM) |



structural formula of EDA-ATP - 5 FAM

Cat. No.: NU-808-5FM

Molecular Formula: C₃₄H₃₂N₇O₂₀P₃ (free acid)

Molecular Weight: 951.58 (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!