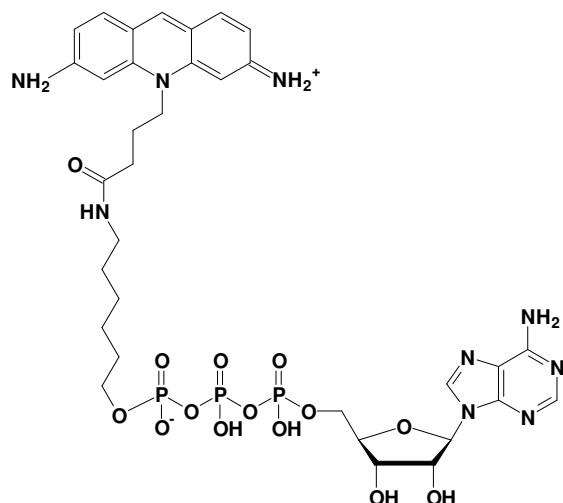


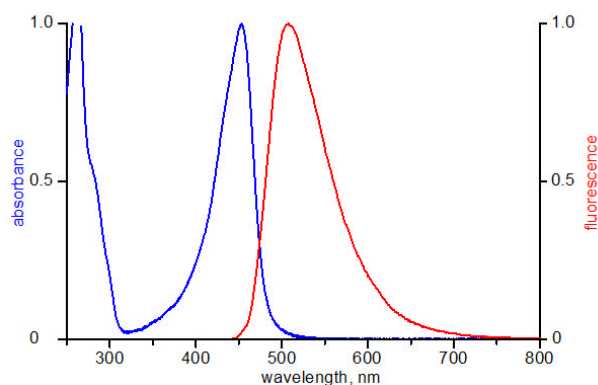
γ -[6-Aminoethyl]-ATP - ATTO 465

γ -[6-Aminoethyl]-adenosine-5'-triphosphate, labeled with ATTO 465, Triethylammonium salt

| Cat. No. | Amount |
|------------|--------------------|
| NU-833-465 | 120 μ l (1 mM) |



structural formula of γ -[6-Aminoethyl]-ATP - ATTO 465



excitation and emission spectrum of ATTO 465

Cat. No.: NU-833-465

Molecular Formula: $C_{33}H_{44}N_9O_{14}P_3$ (free acid)

Molecular Weight: 883.68 (free acid)

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

Spectroscopic properties:

λ_{exc} 453 nm; λ_{em} 508 nm; ϵ 75,000 $cm^{-1} M^{-1}$

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!