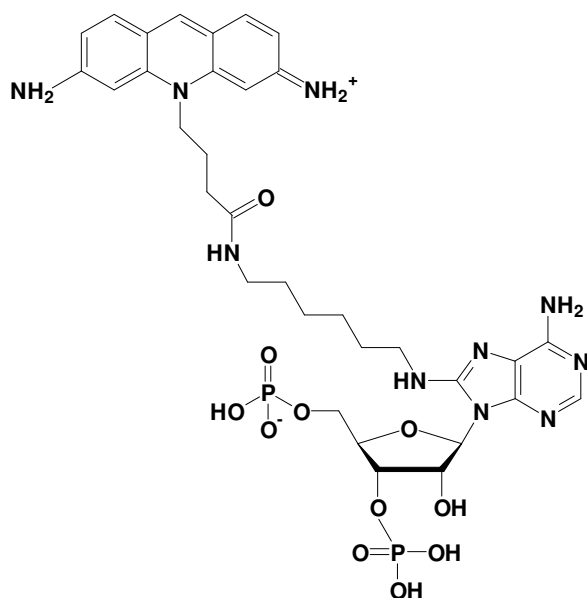


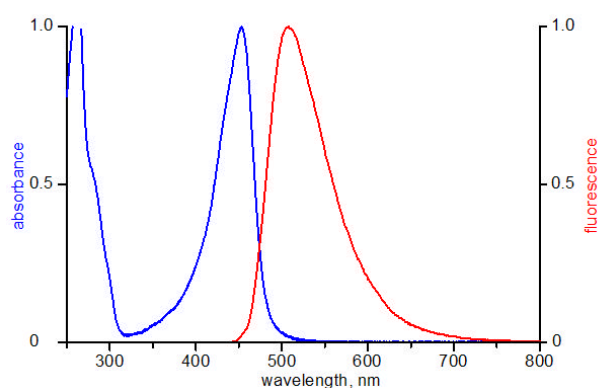
8-[(6-Amino)hexyl]-amino-adenosine-3'/5'-bisphosphate - ATTO 465

8-[(6-Amino)hexyl]-amino-adenosine-3'/5'-bisphosphate, labeled with ATTO 465,
 Triethylammonium salt

Cat. No.	Amount
NU-811-465	120 μ l (1 mM)



structural formula of 8-[(6-Amino)hexyl]-amino-adenosine-3',5'-bisphosphate - ATTO 465



excitation and emission spectrum of ATTO 465

Cat. No.: NU-811-465

Molecular Formula: $C_{33}H_{44}N_{10}O_{11}P_2$ (free acid)

Molecular Weight: 818.72 (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!