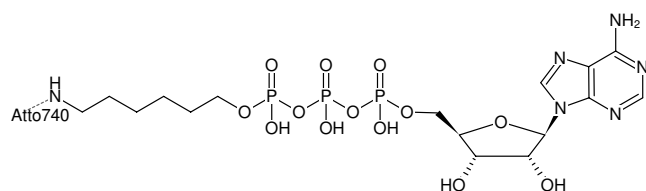




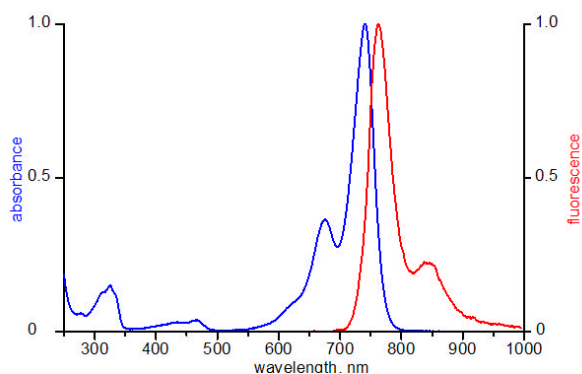
γ -(6-Aminoethyl)-ATP-ATTO-740

γ -(6-Aminoethyl)-adenosine-5'-triphosphate, labeled with ATTO 740, Triethylammonium salt

Cat. No.	Amount
NU-833-740	80 μ l (1 mM)



Structural formula of γ -(6-Aminoethyl)-ATP-ATTO-740



excitation and emission spectrum of ATTO 740

For general laboratory use.

Please centrifuge briefly before opening (volume \leq 2 ml).

Shipping: shipped on gel packs

Storage Conditions: store at -20 °C

Additional Storage Conditions: long time storage at pH 7.5

Short term exposure (up to 1 week cumulative) to ambient temperature possible.

Shelf Life: 12 months after date of delivery

Molecular Formula: $C_{16}H_{29}N_6O_{13}P_3$ - ATTO 740 (free acid)

Molecular Weight: 1055.35 g/mol (free acid)

Purity: \geq 95 % (HPLC)

Form: solution in water

Color: green

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 743 nm, λ_{em} 763 nm, ϵ 120.0 L mmol $^{-1}$ cm $^{-1}$ (Tris-HCl pH 7.5)

Please note: The dye decomposes at pH-levels greater 8.0. The recommended pH for a solution of this product is 7.5.