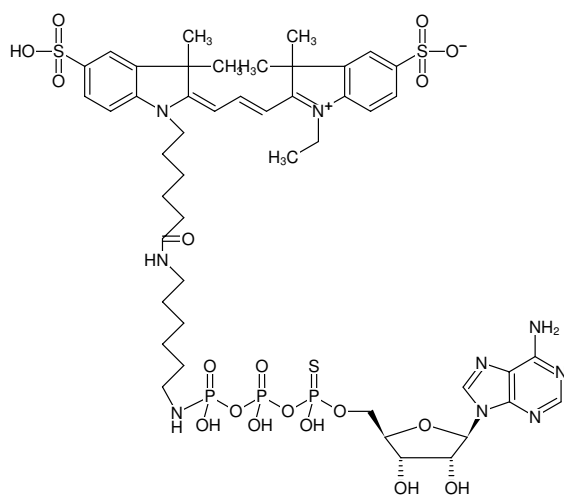




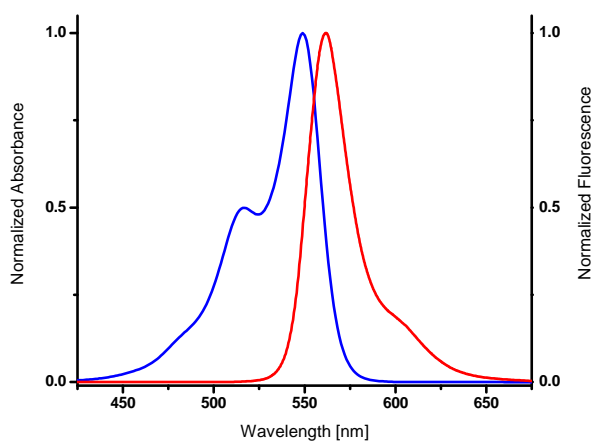
γ -[(6-Aminoethyl)-imido]-ATP α S-Cy3

γ -[(6-Aminoethyl)-imido]-adenosine-5'-(α -thio)-triphosphate, labeled with Cy3, Triethylammonium salt

Cat. No.	Amount
NU-278-CY3	60 μ l (1 mM)



Structural formula of γ -[(6-Aminoethyl)-imido]-ATP α S-Cy3



excitation and emission spectrum of Cy3

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: C₄₇H₆₆N₉O₁₈P₃S₃ (free acid)

Molecular Weight: 1234.19 g/mol (free acid)

Exact Mass: 1233.29 g/mol (free acid)

Purity: \geq 95 % (HPLC)

Form: solution in 10 mM Tris-HCl

Color: pink

Concentration: 1.0 mM - 1.1 mM

pH: 8.0 \pm 0.5

Spectroscopic Properties: λ_{exc} 550 nm, λ_{em} 570 nm, ϵ 150.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)

Please note: This compound contains a phosphoramidite linkage which is hydrolyzed at pH < 7.0.