



EDA-GTP-Cy5

2'/3'-O-(2-Aminoethyl-carbamoyl)-Guanosine-5'-triphosphate, labeled with Cy5, Triethylammonium salt

Cat. No.	Amount
NU-820-CY5	160 µl (1 mM)

For research use only!

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: C₄₆H₆₀N₉O₂₂P₃S₂

Molecular Weight: 1248.07 g/mol (free acid)

Exact Mass: 1247.25 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: solution in water

Color: blue

Concentration: 1.0 mM - 1.1 mM

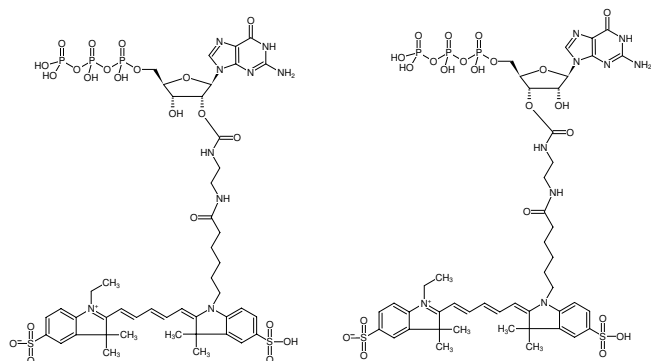
pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 649 nm, λ_{em} 670 nm, ε 250.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)

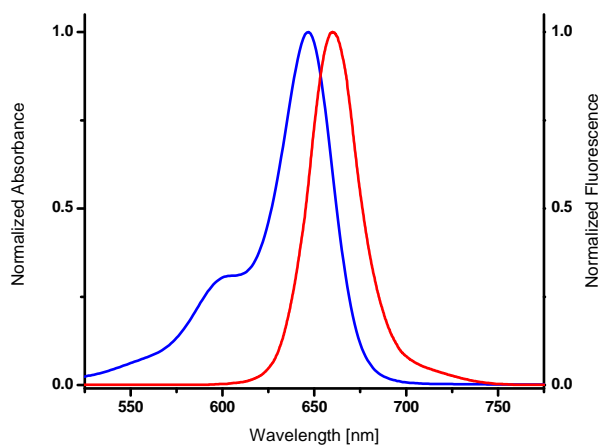
Selected References:

Lang *et al.* (2017) Actin ADP-ribosylation at Threonine148 by Photorhabdus luminescens toxin TccC3 induces aggregation of intracellular F-actin. *Cell. Microbiol.* **19** (1): e12636.

Anderson *et al.* (2011) A fluorescent GTP analog as a specific, high-precision label of microtubules. *BioTechniques* **51** (1):43.



Structural formula of EDA-GTP-Cy5



excitation and emission spectrum of Cy5