

**EDA-m⁷GTP-5-FAM**

2'/3'-O-(2-Aminoethyl-carbamoyl)-7-methyl-guanosine-5'-triphosphate, labeled with 5 FAM, Triethylammonium salt

Cat. No.	Amount
NU-824-5FM	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: C₃₅H₃₄N₇O₂₁P₃ (free acid)

Molecular Weight: 981.61 g/mol (free acid)

Exact Mass: 981.10 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: solution in water

Color: yellow

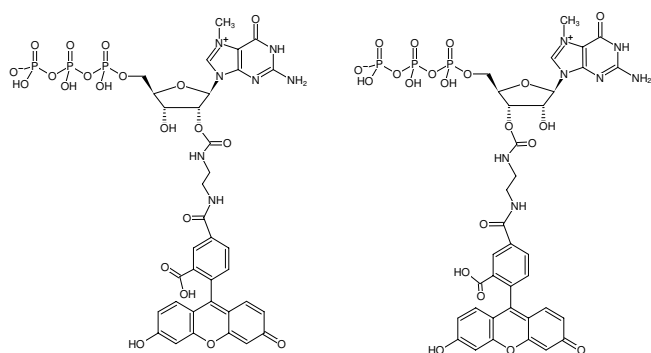
Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

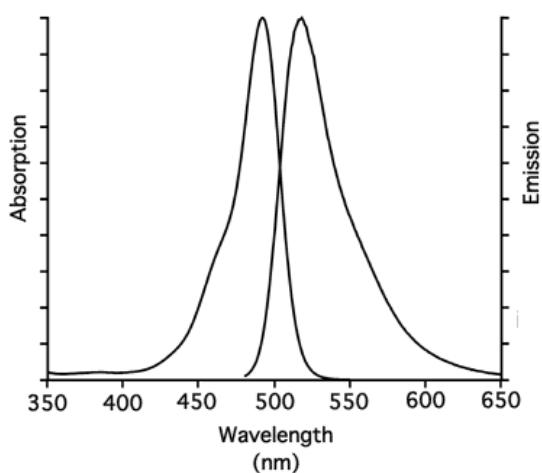
Spectroscopic Properties: λ_{exc} 492 nm, λ_{em} 517 nm, ε 83.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)

Selected References:

Roch *et al.* (2015) An RNA Hybridization Assay for Screening Influenza A Virus Polymerase Inhibitors Using the Entire Ribonucleoprotein Complex. *Assay Drug Dev. Technol.* **13** (8):488.



Structural formula of EDA-m⁷GTP-5-FAM



excitation and emission spectrum of 5-FAM