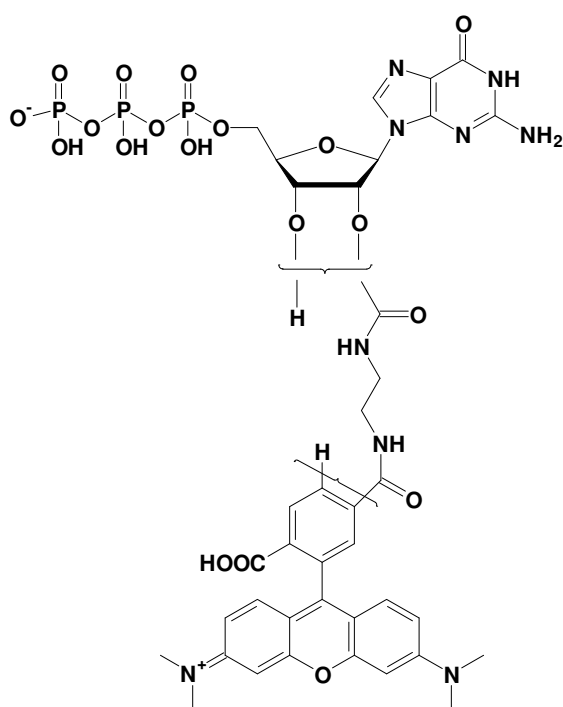


EDA-GTP - 5/6-TAMRA

2'/3'-O-(2-Aminoethyl-carbamoyl)-Guanosine-5'-triphosphate, labeled with 5/6-TAMRA,
 Triethylammonium salt

Cat. No.	Amount
NU-820-TAM	40 μ l / 1 mM



structural formula of EDA-GTP - 5/6-TAMRA

Cat. No.: NU-820-TAM

Molecular Formula: C₃₈H₄₂N₉O₁₉P₃ (free acid)

Molecular Weight: 1021.72 (free acid)

Purity: > 95%, clear aqueous solution, pH 7.5

Spectroscopic properties:

λ_{exc} 545 nm; λ_{em} 575 nm; ϵ 90,000 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!

Selected References:

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

Ahmadian *et. al.* (2005) Monitoring the real-time kinetics of the hydrolysis reaction of guanine nucleotide-binding proteins. *Biol. Chem.* **386**:1105.