



EDA-m⁷GDP-ATTO-465

2'/3'-O-(2-Aminoethyl-carbamoyl)-7-methyl-guanosine-5'-diphosphate, labeled with ATTO 465

Cat. No.	Amount
NU-827-465	120 µl (1 mM)

For research use only!

Shipping: shipped on blue ice

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: C₃₁H₃₈N₁₀O₁₃P₂ (free acid)

Molecular Weight: 820.65 g/mol (free acid)

Exact Mass: 820.21 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: solution in water

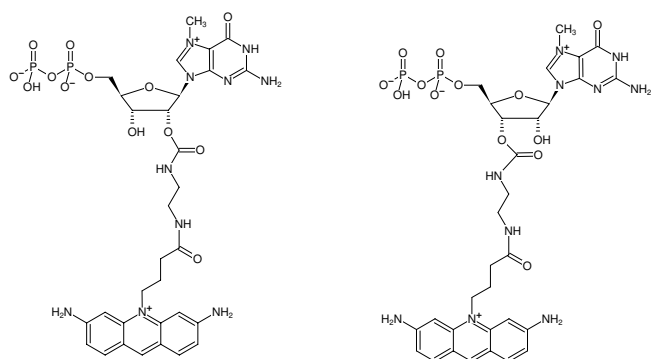
Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

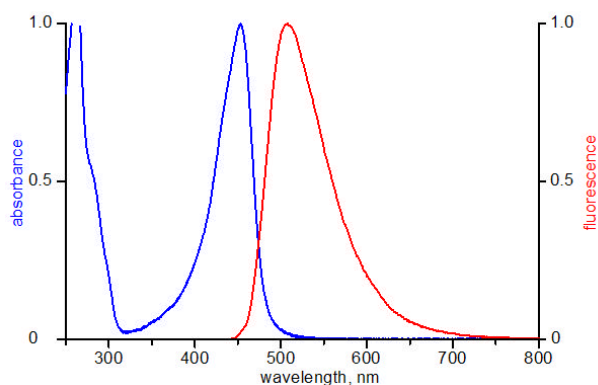
Spectroscopic Properties: λ_{exc} 453 nm, λ_{em} 506 nm, ε 75.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)

Selected References:

Soukarieh *et al.* (2016) Design of nucleotide-mimetic and non-nucleotide inhibitors of the translation initiation factor eIF4E: Synthesis, structural and functional characterisation. *Eur. J. Med. Chem.* **124**:200.



Structural formula of EDA-m⁷GDP-ATTO-465



excitation and emission spectrum of ATTO 465