

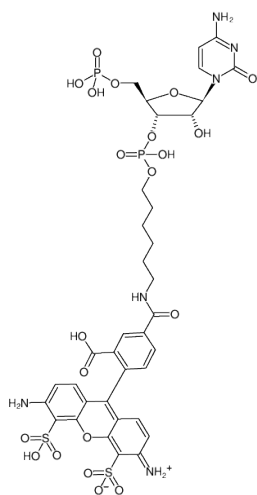


pCp-AZDye488

structural analog to Alexa Fluor® 488

Cytidine-5'-phosphate-3'-(6-aminohexyl)phosphate, labeled with AZDye 488, Triethylammonium salt

Cat. No.	Amount
NU-1706-AZ488	20 µl (1 mM)



Structural formula of pCp-AZDye488

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: C₃₆H₄₀N₆O₂₁P₂S₂ (free acid)

Molecular Weight: 1018.8 g/mol (free acid)

Exact Mass: 1018.12 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

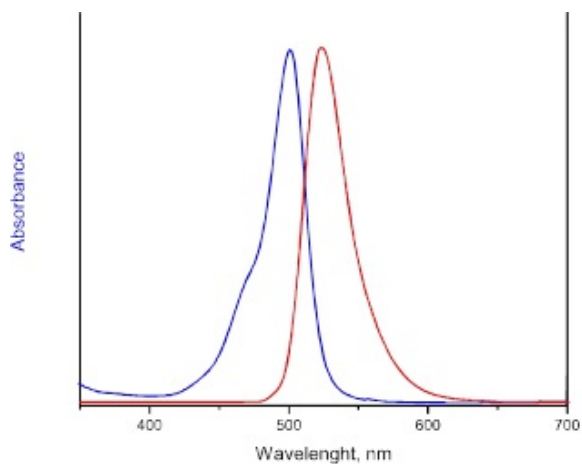
Form: solution in 10 mM Tris-HCl

Color: yellow to orange

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 494 nm, λ_{em} 515 nm, ε 73.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



excitation and emission spectrum of AZDye 488