



zeta-(6-Aminoethyl)-dG6P-AZDye568

structural analog to Alexa Fluor® 568

zeta-(6-Aminoethyl)-2'-deoxyguanosine-5'-hexaphosphate, labeled with AZDye568, Triethylammonium salt

Cat. No.	Amount
NU-284-AZ568-S	20 µl (1 mM)
NU-284-AZ568-L	5 x 20 µl (1 mM)

For general laboratory use.

Please centrifuge briefly before opening (volume ≤2 ml).

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: C₄₉H₆₀N₈O₃₂P₆S₂ (free acid)

Molecular Weight: 1523.01 g/mol (free acid)

Exact Mass: 1522.12 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

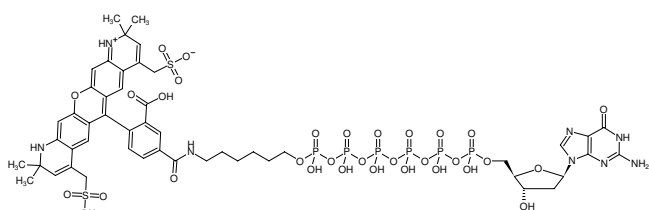
Form: solution in water

Color: red-violet

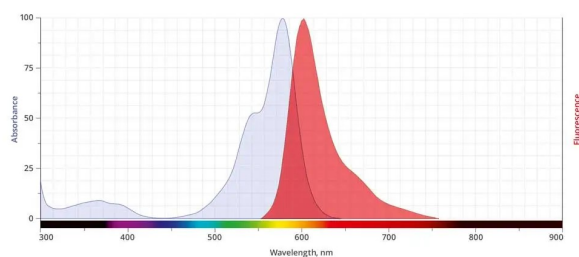
Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ±0.5

Spectroscopic Properties: λ_{exc} 578 nm, λ_{em} 602 nm, ε 88.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



Structural formula of zeta-(6-Aminoethyl)-dG6P-AZDye568



excitation and emission spectrum of AZDye 568