



zeta-(6-Aminohexyl)-dG6P-AZDye647A

structural analog to Alexa Fluor® 647

zeta-(6-Aminohexyl)-2'-deoxyguanosine-5'-hexaphosphate, labeled with AZDye647A, Triethylammonium salt

Cat. No.	Amount
NU-284-AZ647A-S	20 µl (1 mM)
NU-284-AZ647A-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: C₅₂H₇₆N₈O₃₅P₆S₄ (free acid)

Molecular Weight: 1687.28 g/mol (free acid)

Exact Mass: 1686.17 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

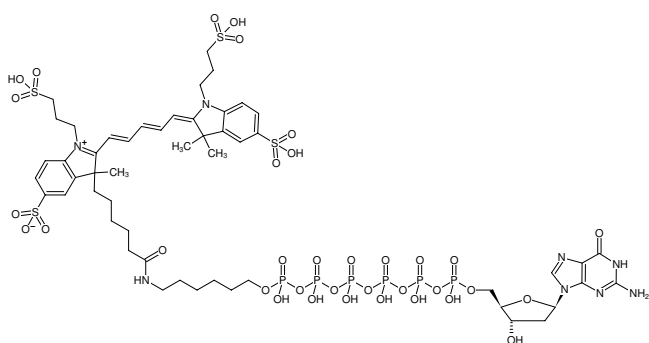
Form: solution in water

Color: blue

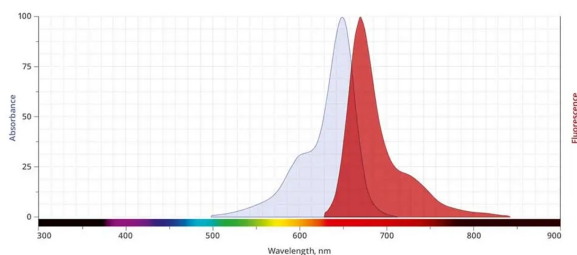
Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ±0.5

Spectroscopic Properties: λ_{exc} 649 nm, λ_{em} 671 nm, ε 270.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



Structural formula of zeta-(6-Aminohexyl)-dG6P-AZDye647A



excitation and emission spectrum of AZDye 647A