



## zeta-(6-Aminohexyl)-dT6P-AZDye555A

structural analog to Alexa Fluor® 555

zeta-(6-Aminohexyl)-2'-deoxythymidine-5'-hexaphosphate, labeled with AZDye555A, Triethylammonium salt

Cat. No.	Amount
NU-285-AZ555A-S	20 µl (1 mM)
NU-285-AZ555A-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:** C<sub>50</sub>H<sub>75</sub>N<sub>5</sub>O<sub>36</sub>P<sub>6</sub>S<sub>4</sub> (free acid)

**Molecular Weight:** 1636.23 g/mol (free acid)

**Exact Mass:** 1635.15 g/mol (free acid)

**Purity:** ≥ 95 % (HPLC)

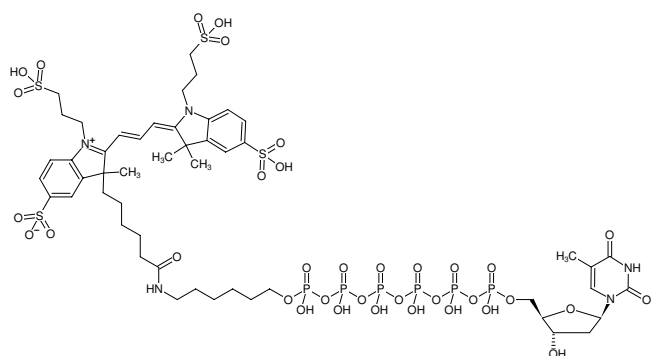
**Form:** solution in water

**Color:** red

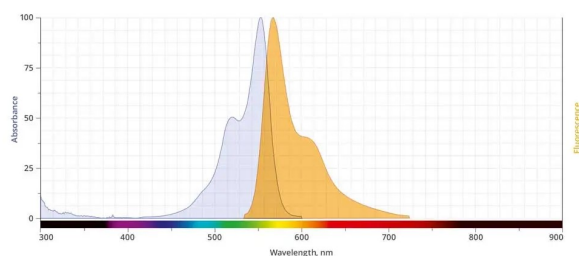
**Concentration:** 1.0 mM - 1.1 mM

**pH:** 7.5 ±0.5

**Spectroscopic Properties:** λ<sub>exc</sub> 555 nm, λ<sub>em</sub> 572 nm, ε 155.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)



Structural formula of zeta-(6-Aminohexyl)-dT6P-AZDye555A



excitation and emission spectrum of AZDye 555A