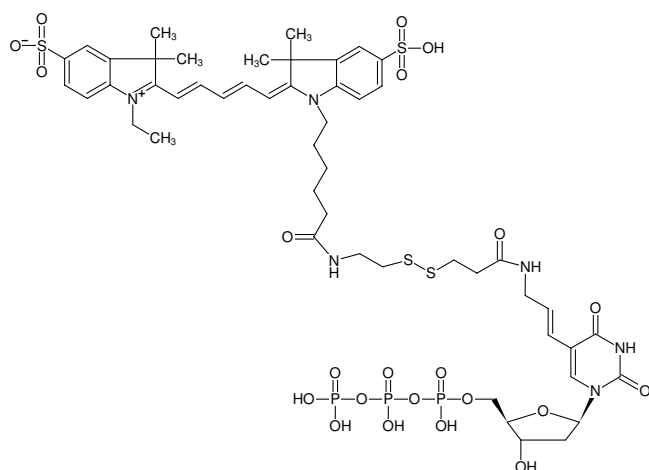




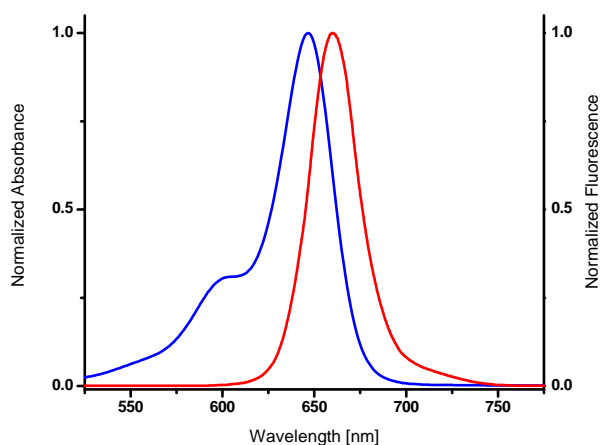
5-Aminoallyl-dUTP-SS-Cy5

dUTP-SS-Cy5
Triethylammonium salt

Cat. No.	Amount
NU-803-SS-CY5	20 µl (1 mM)



Structural formula of 5-Aminoallyl-dUTP-SS-Cy5



excitation and emission spectrum of Cy5

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: C₅₀H₆₇N₆O₂₂P₃S₄ (free acid)

Molecular Weight: 1325.27 g/mol (free acid)

Exact Mass: 1324.24 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} 670 nm, ε 250.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)

Selected References:

Goltsev *et al.* (2018) Deep Profiling of Mouse Splenic Architecture with CODEX Multiplexed Imaging. *Cell*. **174** (4):968-981.e15.