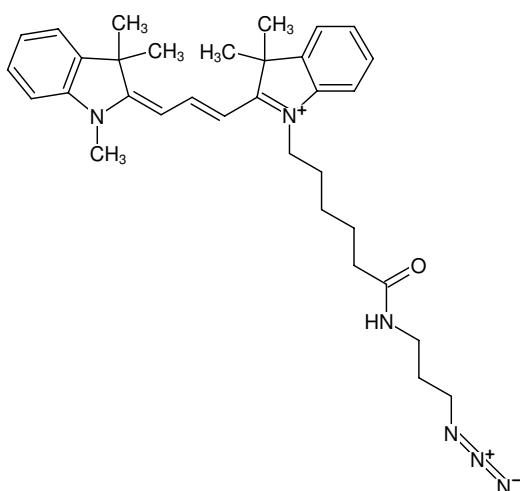




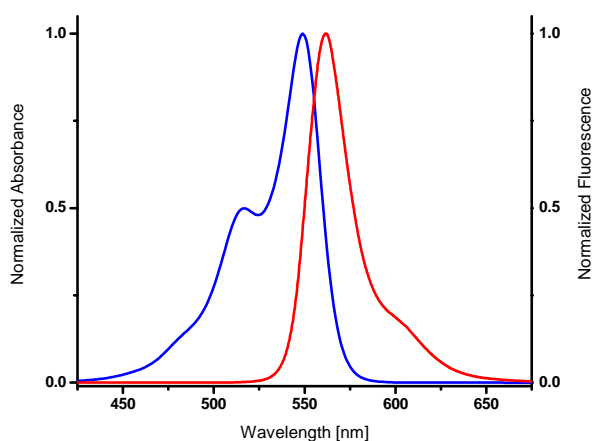
Cy3-Azide

Abs/Em = 555/570 nm

Cat. No.	Amount
CLK-046-1	1 mg
CLK-046-5	5 x 1 mg



Structural formula of Cy3-Azide



excitation and emission spectrum of Cy3

For general laboratory use.

Shipping: shipped at ambient temperature

Storage Conditions: store at -20 °C

Additional Storage Conditions: store dark and dry

Shelf Life: 12 months after date of delivery

Molecular Formula: C₃₃H₄₃N₆O (free cation)

Molecular Weight: 539.75 g/mol (free cation)

Exact Mass: 539.35 g/mol (free cation)

Purity: ≥ 90 % (HPLC)

Form: powder

Color: red

Solubility: soluble in DMF, DMSO, Dichlormethan
insoluble in water

Spectroscopic Properties: λ_{abs} 555 nm, λ_{em} 570 nm,
ε 150.0 L mmol⁻¹ cm⁻¹ (pH 9)