



DBCO-Sulfo-Cy5

Abs/Em = 646/661 nm

Dibenzylcyclooctyne-Sulfo-Cy5

Cat. No.	Amount
CLK-A130-1	1 mg
CLK-A130-5	5 x 1 mg
CLK-A130-25	25 mg

For general laboratory use.

Shipping: shipped at ambient temperature

Storage Conditions: store at -20 °C

Shelf Life: 12 months after date of delivery

Molecular Formula: C₅₂H₅₆N₄O₁₁S₃ (free acid)

Molecular Weight: 1009.22 g/mol (free acid)

Exact Mass: 1008.31 g/mol (free acid)

CAS#: 1564286-24-3 (inner salt)

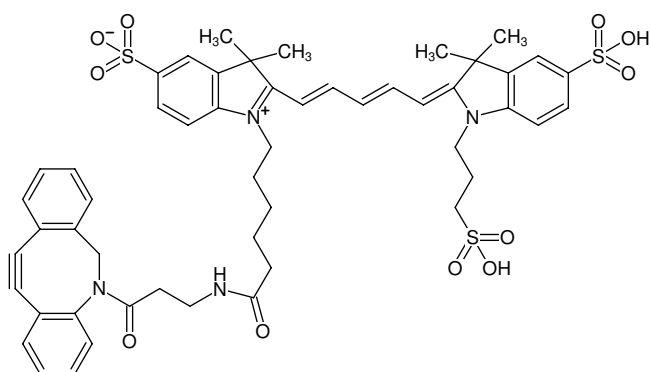
Purity: ≥ 90 % (HPLC)

Form: powder

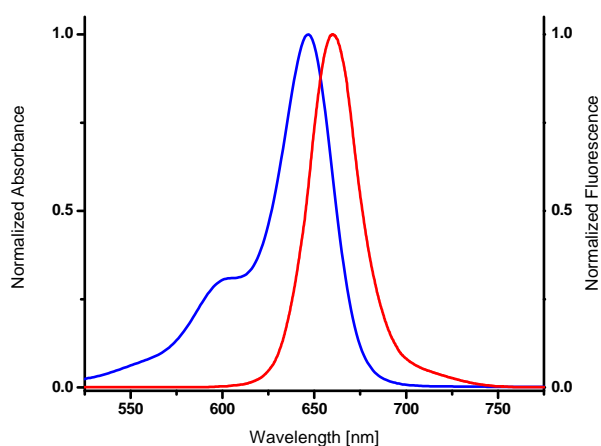
Color: blue

Solubility: DMF, DMSO, water

Spectroscopic Properties: λ_{abs} 646 nm, λ_{em} 661 nm, ε 251.0 L mmol⁻¹ cm⁻¹



Structural formula of DBCO-Sulfo-Cy5



excitation and emission spectrum of Cy5