



DBCO-Sulfo-Cy3

Abs/Em = 553/563 nm
Dibenzylcyclooctyne-Sulfo-Cy3

Cat. No.	Amount
CLK-A140-1	1 mg
CLK-A140-5	5 x 1 mg
CLK-A140-25	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: C₅₀H₅₄N₄O₁₁S₃ (free acid)

Molecular Weight: 983.18 g/mol (free acid)

Exact Mass: 982.30 g/mol (free acid)

CAS#: 1782950-79-1

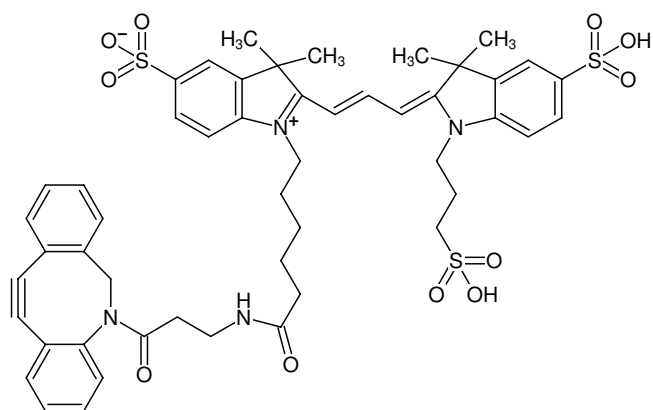
Purity: ≥ 90 % (HPLC)

Form: solid

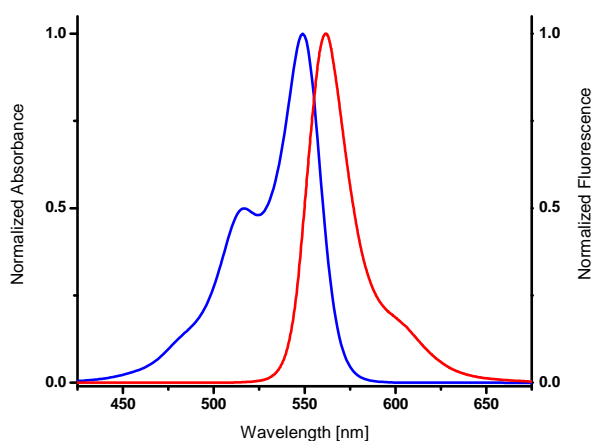
Color: red

Solubility: DMF, DMSO, water

Spectroscopic Properties: λ_{abs} 553 nm, λ_{em} 563 nm, ε 151.0 L mmol⁻¹ cm⁻¹



Structural formula of DBCO-Sulfo-Cy3



excitation and emission spectrum of Cy3