



DBCO-PEG₄-5/6-Carboxyrhodamine 110

Abs/Em = 501/526 nm
 Dibenzylcyclooctyne-PEG₄-5/6-Carboxyrhodamine 110
 DBCO-PEG₄-Rhodamine Green

Cat. No.	Amount
CLK-A127-1	1 mg
CLK-A127-5	5 x 1 mg
CLK-A127-25	25 mg

For research use only!

Shipping: shipped at ambient temperature

Storage Conditions: store at -20 °C

Shelf Life: 12 months after date of delivery

Molecular Formula: C₅₀H₄₉N₅O₁₀

Molecular Weight: 879.95 g/mol

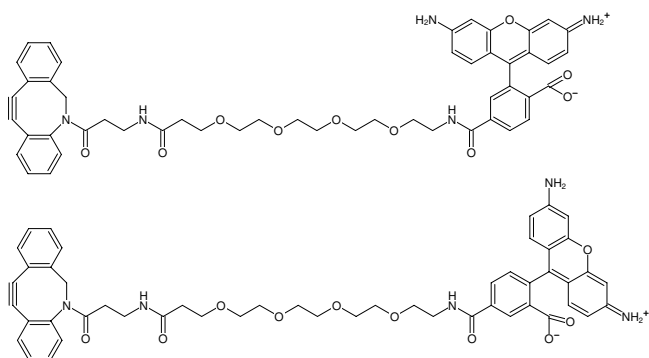
Purity: > 90 % (HPLC)

Form: solid

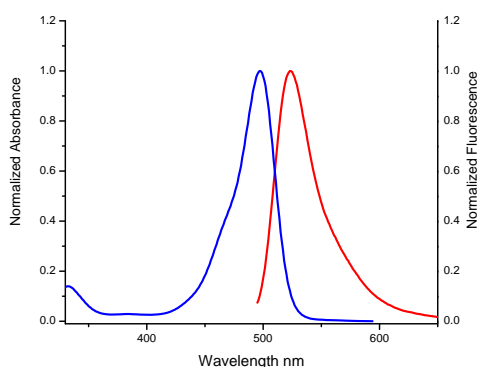
Color: red

Solubility: DMF, DMSO

Spectroscopic Properties: λ_{abs} 501 nm, λ_{em} 526 nm,
 ε 74.0 L mmol⁻¹ cm⁻¹



Structural formula of DBCO-PEG₄-5/6-Carboxyrhodamine 110



excitation and emission spectrum of 5/6-Carboxyrhodamine 110