



DBCO-AF488

Abs/Em = 494/517 nm
also known as Alexa Fluor® 488-DBCO

Cat. No.	Amount
CLK-1278-1	1 mg
CLK-1278-5	5 x 1 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: C₃₉H₂₈N₄O₁₁S₂

Molecular Weight: 792.79 g/mol

Exact Mass: 792.12 g/mol

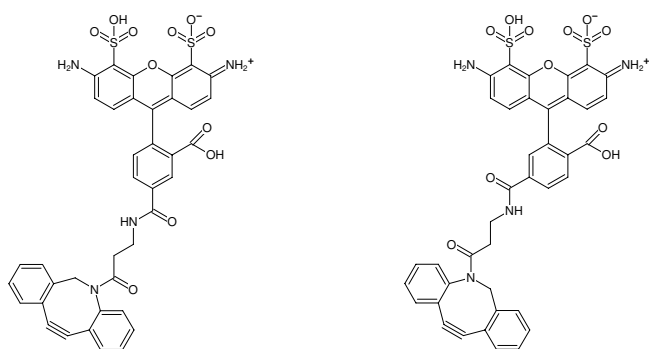
Purity: > 95 % (HPLC)

Form: solid

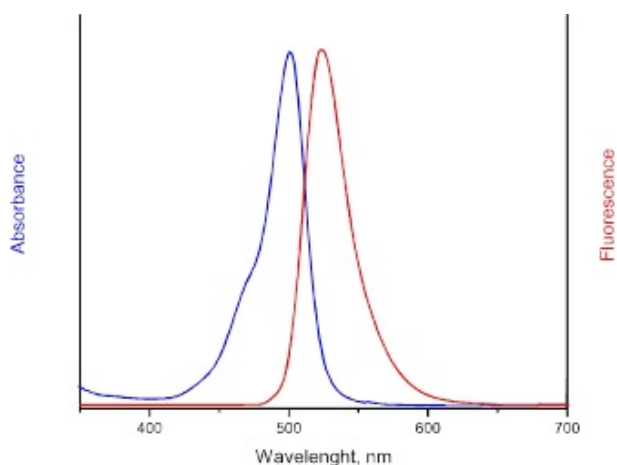
Color: orange to red

Solubility: DMSO, DMF

Spectroscopic Properties: λ_{exc} 494 nm, λ_{em} 517 nm, ε 73.0 L mmol⁻¹ cm⁻¹



Structural formula of DBCO-AF488



excitation and emission spectrum of AF488