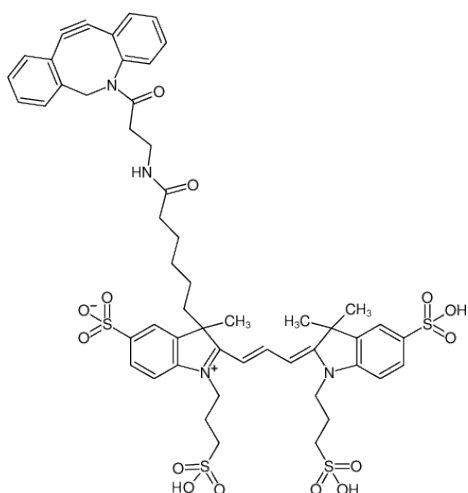


**DBCO-AF555**

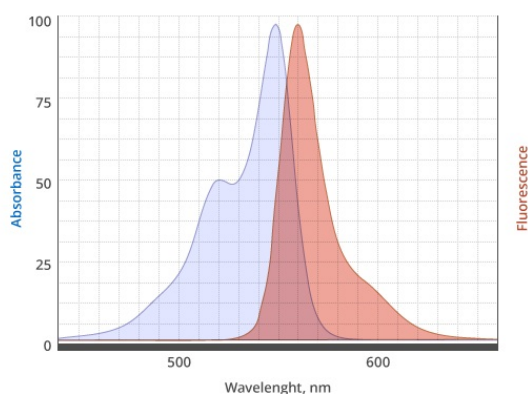
Abs/Em = 555/572 nm

structural analog to Alexa Fluor® 555

Cat. No.	Amount
CLK-1290-1	1 mg
CLK-1290-5	5 x 1 mg



Structural formula of DBCO-AF555

**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:** C<sub>52</sub>H<sub>58</sub>N<sub>4</sub>O<sub>14</sub>S<sub>4</sub>**Molecular Weight:** 1091.29 g/mol**Exact Mass:** 1090.28 g/mol**Purity:** ≥ 90 % (HPLC)**Form:** solid**Color:** red**Solubility:** DMSO, DMF**Spectroscopic Properties:** λ<sub>exc</sub> 555 nm, λ<sub>em</sub> 572 nm, ε 155.0 L mmol<sup>-1</sup> cm<sup>-1</sup>

excitation and emission spectrum of AF555