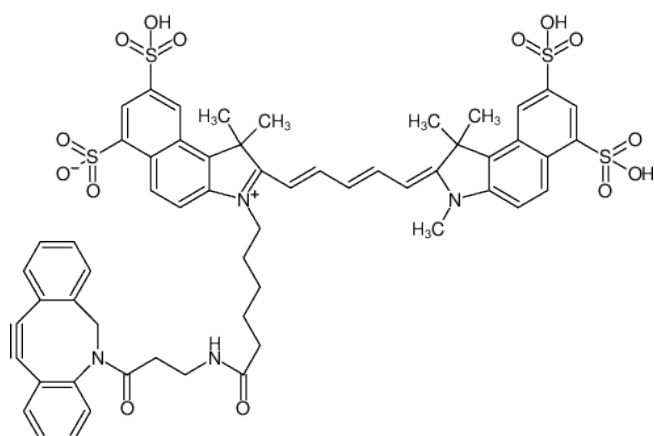


**DBCO-Cy5.5**

Abs/Em = 678/694 nm  
 Dibenzylcyclooctyne-Cy5.5  
 Triethylammonium salt

Cat. No.	Amount
CLK-1046A-1	1 mg
CLK-1046A-5	5 x 1 mg
CLK-1046A-25	25 mg



Structural formula of DBCO-Cy5.5

**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>58</sub>H<sub>56</sub>N<sub>4</sub>O<sub>14</sub>S<sub>4</sub>**Molecular Weight:** 1161.34 g/mol**Exact Mass:** 1160.26 g/mol**Purity:** ≥ 90 % (HPLC)**Form:** solid**Color:** blue**Solubility:** DMF, DMSO, water**Spectroscopic Properties:** λ<sub>abs</sub> 678 nm, λ<sub>em</sub> 694 nm,  
ε 190.0 L mmol<sup>-1</sup> cm<sup>-1</sup>