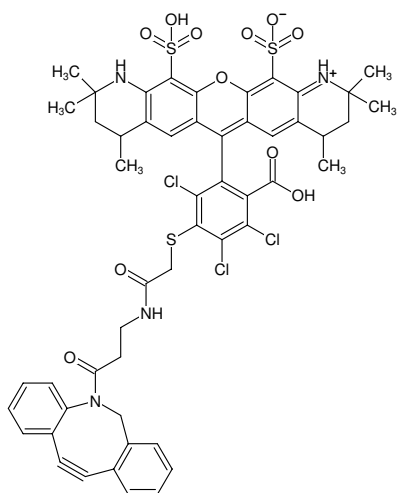




## DBCO-AF546

Abs/Em = 554/570 nm  
also known as Alexa Fluor® 546-DBCO

Cat. No.	Amount
CLK-1286-1	1 mg
CLK-1286-5	5 x 1 mg



Structural formula of DBCO-AF546

**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>47</sub>Cl<sub>3</sub>N<sub>4</sub>O<sub>11</sub>S<sub>3</sub>

**Molecular Weight:** 1106.50 g/mol

**Exact Mass:** 1104.15 g/mol

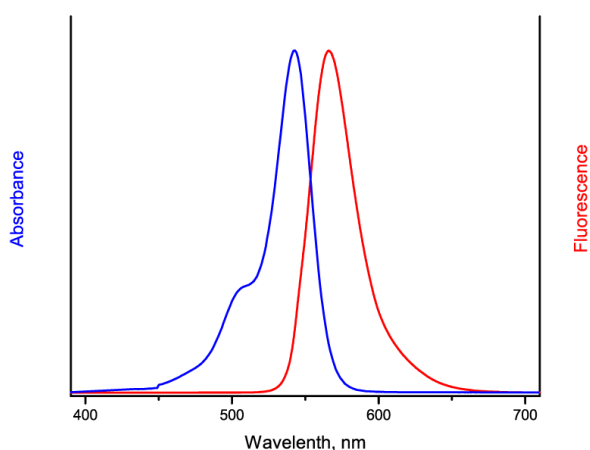
**Purity:** ≥ 90 % (HPLC)

**Form:** solid

**Color:** red

**Solubility:** DMSO, DMF

**Spectroscopic Properties:** λ<sub>exc</sub> 554 nm, λ<sub>em</sub> 570 nm, ε 110.0 L mmol<sup>-1</sup> cm<sup>-1</sup>



excitation and emission spectrum of AF546