



## DBCO-AF594

Abs/Em = 590/617 nm  
also known as Alexa Fluor® 594-DBCO

Cat. No.	Amount
CLK-1298-1	1 mg
CLK-1298-5	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:** C<sub>53</sub>H<sub>48</sub>N<sub>4</sub>O<sub>11</sub>S<sub>2</sub>

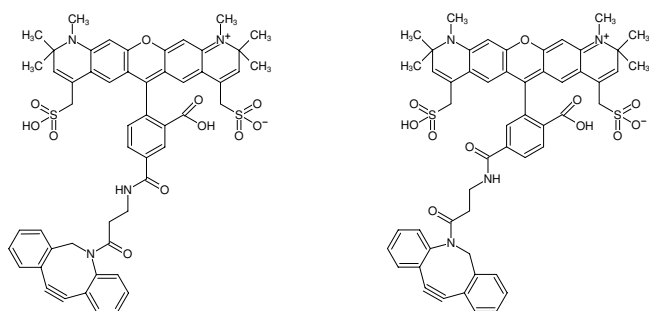
**Molecular Weight:** 981.10 g/mol

**Purity:** > 95 % (HPLC)

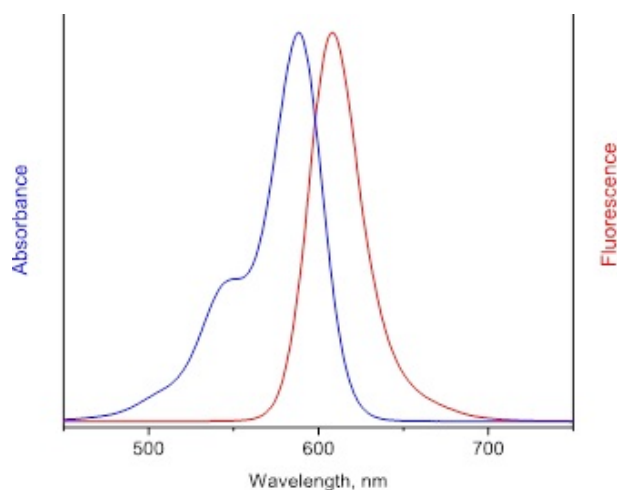
**Form:** blue solid

**Solubility:** DMSO, DMF

**Spectroscopic Properties:** λ<sub>exc</sub> 590 nm, λ<sub>em</sub> 617 nm, ε 92.0 L mmol<sup>-1</sup> cm<sup>-1</sup>



Structural formula of DBCO-AF594



excitation and emission spectrum of AF594