



## 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:** C<sub>39</sub>H<sub>28</sub>N<sub>4</sub>O<sub>11</sub>S<sub>2</sub>

**Molecular Weight:** 792.79 g/mol

**Exact Mass:** 792.12 g/mol

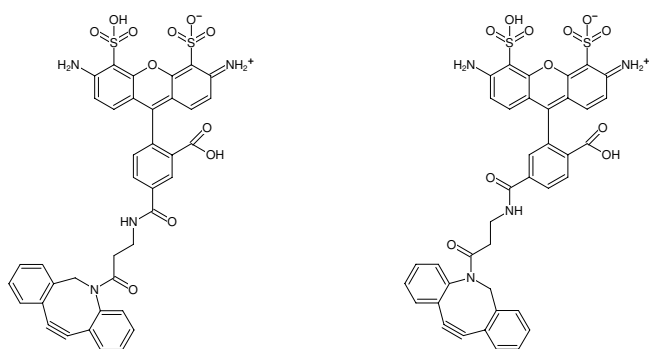
**Purity:** > 95 % (HPLC)

**Form:** solid

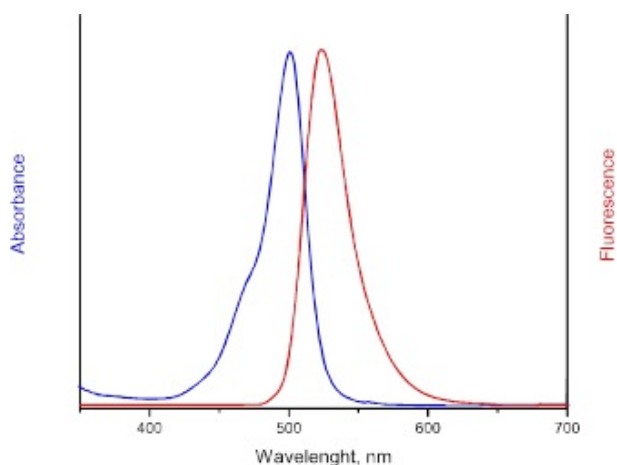
**Color:** orange to red

**Solubility:** DMSO, DMF

**Spectroscopic Properties:** λ<sub>exc</sub> 494 nm, λ<sub>em</sub> 517 nm, ε 73.0 L mmol<sup>-1</sup> cm<sup>-1</sup>



Structural formula of DBCO-AF488



excitation and emission spectrum of AF488