



## AF647-Picolyl-Azide

Abs/Em = 648/671 nm  
structural analog to Alexa Fluor® 647

Cat. No.	Amount
CLK-1300-1	1 mg
CLK-1300-5	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>46</sub>H<sub>58</sub>N<sub>8</sub>O<sub>14</sub>S<sub>4</sub>

**Molecular Weight:** 1077.27 g/mol

**Exact Mass:** 1076.31 g/mol

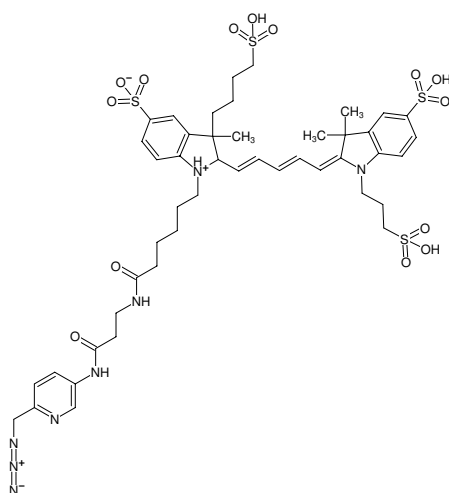
**Purity:** > 95 % (HPLC)

**Form:** solid

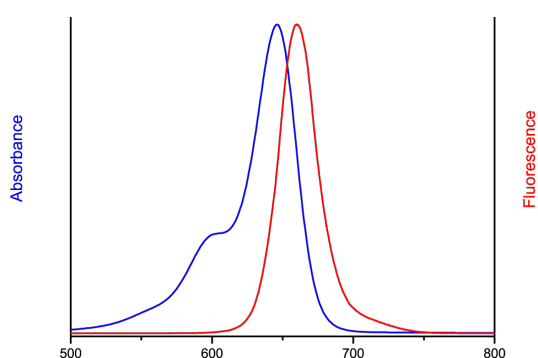
**Color:** blue

**Solubility:** DMSO, DMF

**Spectroscopic Properties:** λ<sub>exc</sub> 648 nm, λ<sub>em</sub> 671 nm, ε 270.0 L mmol<sup>-1</sup> cm<sup>-1</sup>



Structural formula of AF647-Picolyl-Azide



excitation and emission spectrum of AF647