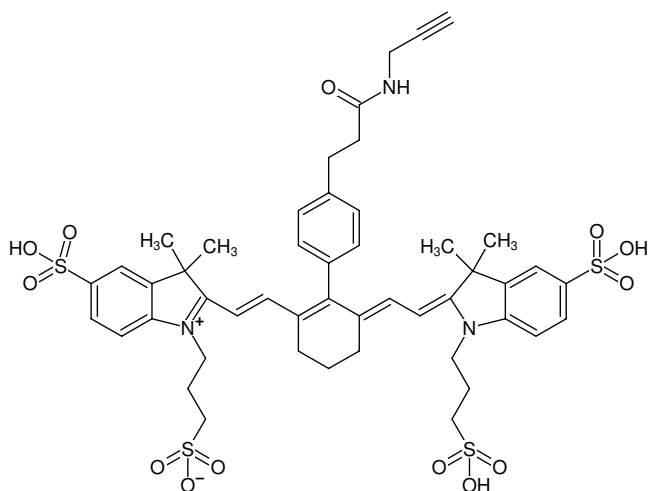




## Cy7-Alkyne

Abs/Em = 753/775 nm  
Triethylammonium salt

Cat. No.	Amount
CLK-1053-1	1 mg
CLK-1053-5	5 x 1 mg
CLK-1053-25	25 mg



Structural formula of Cy7-Alkyne

**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>48</sub>H<sub>55</sub>N<sub>3</sub>O<sub>13</sub>S<sub>4</sub>

**Molecular Weight:** 1010.22 g/mol

**Purity:** > 90 % (HPLC)

**Form:** solid

**Color:** green

**Solubility:** DMF, DMSO, MeOH, water

**Spectroscopic Properties:** λ<sub>abs</sub> 753 nm, λ<sub>em</sub> 775 nm,  
ε 255.0 L mmol<sup>-1</sup> cm<sup>-1</sup>