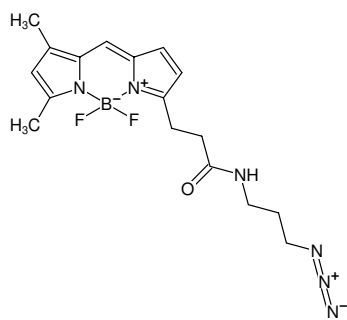


**Azide-BDP-FL**

Abs/Em = 503/512 nm  
also known as Azide-BODIPY® FL

Cat. No.	Amount
CLK-044-1	1 mg
CLK-044-5	5 x 1 mg



Structural formula of Azide-BDP-FL

**For research use only!**

**Shipping:** shipped at ambient temperature

**Storage Conditions:** store at -20 °C

**Additional Storage Conditions:** store dark and dry

**Shelf Life:** 12 months after date of delivery

**Molecular Formula:** C<sub>17</sub>H<sub>21</sub>BF<sub>2</sub>N<sub>6</sub>O

**Molecular Weight:** 374.18 g/mol

**Purity:** > 95 % (HPLC-MS, H NMR)

**Form:** orange solid

**Solubility:** alcohols, DMF, DMSO

**Spectroscopic Properties:** λ<sub>exc</sub> 503 nm, λ<sub>em</sub> 512 nm,  
ε 80.0 L mmol<sup>-1</sup> cm<sup>-1</sup>