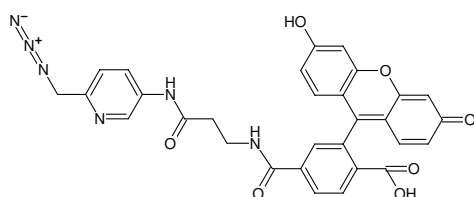
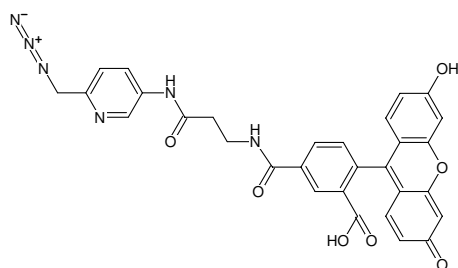




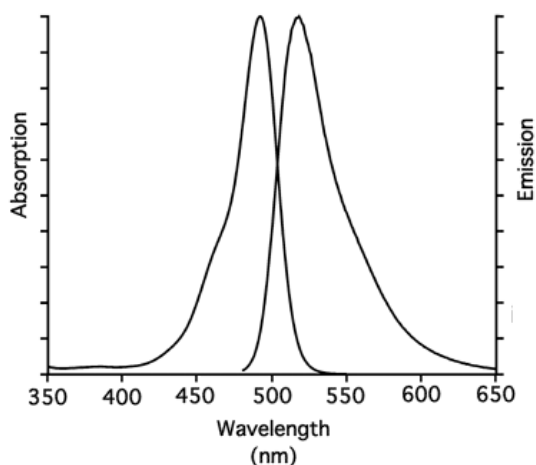
### Picolyl-Azide-5/6-FAM

Abs/Em = 490/510 nm

Cat. No.	Amount
CLK-1180-1	1 mg
CLK-1180-5	5x 1 mg



Structural formula of Picolyl-Azide-5/6-FAM



excitation and emission spectrum of 5/6-FAM

**For general laboratory use.**

**Shipping:** shipped at ambient temperature

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

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

**Molecular Formula:** C<sub>30</sub>H<sub>22</sub>N<sub>6</sub>O<sub>7</sub> (free acid)

**Molecular Weight:** 578.53 g / mol (free acid)

**Exact Mass:** 578.15 g / mol (free acid)

**Purity:** ≥ 90 % (HPLC)

**Form:** solid

**Color:** yellow

**Solubility:** DMSO, DMF

**Spectroscopic Properties:** λ<sub>exc</sub> 490 nm, λ<sub>em</sub> 510 nm, ε 76.0 L mmol<sup>-1</sup> cm<sup>-1</sup>

**Selected References:**

Jiang *et al.* (2014) Monitoring Dynamic Glycosylation in Vivo Using Supersensitive Click Chemistry. *Bioconj. Chem.* **25** (4):698.

Uttamapinant *et al.* (2012) Fast, cell-compatible click chemistry with copper-chelating azides for biomolecular labeling. *Angew. Chem. Int. Ed. Engl.* **51** (24):5852.