

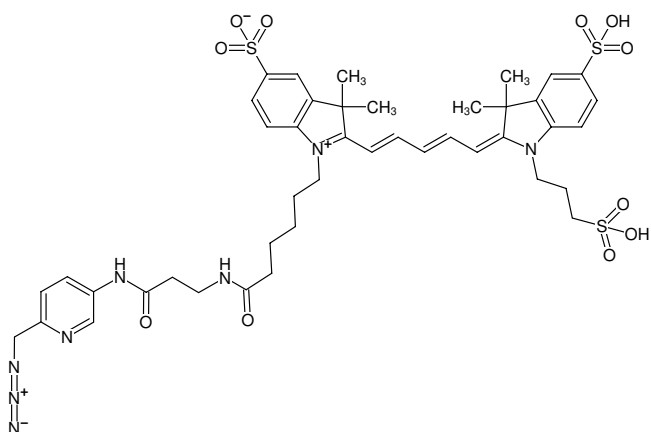

**Picolyl-Azide-Sulfo-Cy5**

Abs/Em = 647/663 nm

Cat. No.	Amount
CLK-1177-1	1 mg
CLK-1177-5	5 x 1 mg
CLK-1177-25	25 mg

**Selected References:**

 Jiang *et al.* (2014) Monitoring Dynamic Glycosylation in Vivo Using Supersensitive Click Chemistry. *Bioconjug. 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.


Structural formula of Picolyl-Azide-Sulfo-Cy5

**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>43</sub>H<sub>52</sub>N<sub>8</sub>O<sub>11</sub>S<sub>3</sub> (free acid)

**Molecular Weight:** 953.11 g / mol (free acid)

**Exact Mass:** 952.29 g / mol (free acid)

**Purity:** ≥ 90 % (HPLC)

**Form:** solid

**Color:** blue

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

**Spectroscopic Properties:** λ<sub>exc</sub> 647 nm, λ<sub>em</sub> 663 nm, ε 255.0 L mmol<sup>-1</sup> cm<sup>-1</sup>