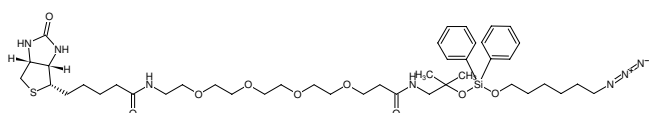


**DADPS Biotin-Azide**

Cat. No.	Amount
CLK-1330-5	5 mg
CLK-1330-25	25 mg



Structural formula of DADPS Biotin-Azide

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:** C₄₃H₆₇N₇O₉Si**Molecular Weight:** 886.19 g/mol**Exact Mass:** 885.45 g/mol**Purity:** ≥ 90 % by reaction with DBCO-compound followed by cleavage**Form:** oil to grey amorphous solid**Solubility:** DMSO, DMF