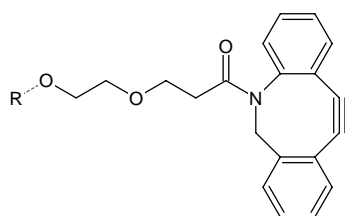




DBCO Magnetic Beads

Cat. No.	Amount
CLK-1037-1	1 ml



Structural formula of DBCO Magnetic Beads

For research use only!**Shipping:** shipped at ambient temperature**Storage Conditions:** store at 4 °C**Additional Storage Conditions:** do not freeze**Shelf Life:** 12 months after date of delivery**Form:** 10 mg/ml suspension containing 0.05 % Kathon**Applications:**

DBCO Magnetic Beads are an efficient matrix to covalently capture CLICK-functionalized proteins by Cu(I)-free click chemistry reaction. The proteins of interest need to be metabolically, enzymatically or chemically azide-tagged.

Subsequently, the DBCO Magnetic Beads containing the covalently attached proteins can be washed with high stringency, virtually eliminating any non-specifically bound proteins.

DBCO Magnetic Beads are not stable in organic solvents and high urea concentrations.

Degree of substitution: 30-50 nmol DBCO-groups/mg resin**Surface area:** $5.7 \cdot 10^9 \mu\text{m}^2/\text{mg}$ **Mean bead diameter:** 0.97 μm **Density:** 1.33 g/cm³**Polymere composition:** benzene, ethynyl- and homopolymer (encapsulated magnetite)