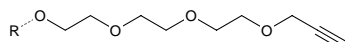




Alkyne Magnetic Beads

Cat. No.	Amount
CLK-1035-1	1 ml



Structural formula of Alkyne 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:

Alkyne Magnetic Beads are an efficient magnetic matrix to covalently capture CLICK-functionalized proteins by a Cu(I)-catalyzed azide-alkyne cycloaddition (CuAAC). The proteins of interest need to be metabolically, enzymatically or chemically azide-tagged. Subsequently, the azide magnetic beads containing the covalently attached proteins can be washed with high stringency, virtually eliminating any non-specifically bound proteins.

Alkyne Magnetic Beads are not stable in organic solvents and high urea concentrations. Use Alkyne Agarose (CLK-1032) for mass spectrometry experiments.

Degree of substitution: 30-50 nmol Alkyne-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/cm^3

Polymere composition: benzene, ethylenyl- and homopolymer (encapsulated magnetite)