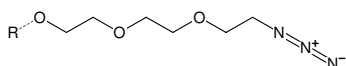




Azide Agarose

Cat. No.	Amount
CLK-1038-2	2 ml



Structural formula of Azide Agarose

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: 50 % aqueous suspension containing 30 % EtOH

Color: off-white

Applications:

Azide Agarose is an efficient matrix to covalently capture CLICK-functionalized proteins by a Cu(I)-catalyzed azide-alkyne cycloaddition (CuAAC) or a Cu(I)-free click chemistry reaction. The proteins of interest need to be metabolically, enzymatically or chemically alkyne-tagged (Cu(I)-catalyzed reaction) or DBCO-tagged (Cu(I)-free reaction).

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

For a comprehensive protocol for Cu(I)-catalyzed purification of alkyne-tagged proteins, please refer to our Click Chemistry Capture Kit (CLK-1065).

Degree of substitution: 10-20 µmol Azide-groups/ml resin

Matrix: 6 % cross-linked agarose

Bead size: 50-150 µm, spherical