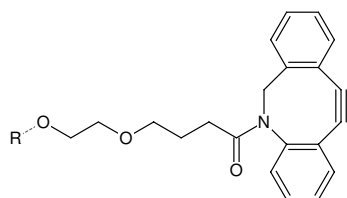


**DBCO Agarose**

Cat. No.	Amount
CLK-1034-2	2 ml



Structural formula of DBCO 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 % slightly yellow to yellow aqueous suspension containing 30 % EtOH**Applications:**

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

Degree of substitution: 10-20 µmol DBCO-groups/ml resin**Matrix:** 6 % cross-linked agarose**Bead size:** 50-150 µm, spherical