



Dde Alkyne Agarose

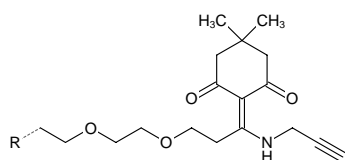
agarose resin with alkyne groups and cleavable Dde linker

Cat. No.	Amount
CLK-1140-2	2 ml

Degree of substitution: 5 - 20 μmol Alkyne-groups/ml resin

Matrix: 6 % cross-linked agarose

Bead size: 50-150 μm ; spherical



Structural formula of Dde Alkyne 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 % off-white aqueous suspension containing 20 % EtOH