



### PC Alkyne Agarose

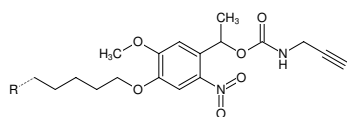
agarose resin with alkyne groups and photocleavable linker

Cat. No.	Amount
CLK-1142-2	2 ml

**Degree of substitution:** 5 - 20  $\mu\text{mol}$  Alkyne-groups/ml resin

**Matrix:** 6 % cross-linked agarose

**Bead size:** 50-150  $\mu\text{m}$ , spherical



Structural formula of PC 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 % aqueous suspension containing 20 % EtOH

**Color:** off-white