

Agarose characteristics

Bead/Particle size

45-165 μm

Recommended linear flow rate

11.5 cm/h

Maximum pressure

0.25 bar (3.6 psi)

pH stability

short term: 4 - 9 / long term: 7.5

Chemical stability

Stable to all solutions commonly used in gel filtration including 8 M urea and 6 M guanidine hydrochloride

Not stable in organic solvents!

Sterilization

Not autoclavable!

R= Agarose