



Dear Extra Fine P-Gel User,

This extra fine P-gel is a specially formulated polyacrylamide material for high resolution gel filtration. The size specification of beads, $< 45 \mu\text{m}$ in diameter, allows higher resolution, with a corresponding decrease in flow rate. To improve flow rates, we recommend removal of the smaller particles following swelling. Gently suspend the gel in a two- to four-fold excess of eluant. Allow 85-95% of the gel to settle out. Remove the remaining gel and supernatant rapidly with suction. Repeat twice, or until a rapidly sedimenting bed is obtained.