

Protein Purification Flowchart Using StrepCaptureXP Resin FF

1 Sample Preparation

Ensure the sample is clear and filtered (0.22 μm) before loading. Use your standard protocol.

Tip: Keep cold. Add protease inhibitors if needed.



-  Resin
-  Target protein
-  Impurities

2 Column Chromatography (Gravity Column)

A. Column Packing & Equilibration: Equilibrate with 10 bed volumes washing buffer.

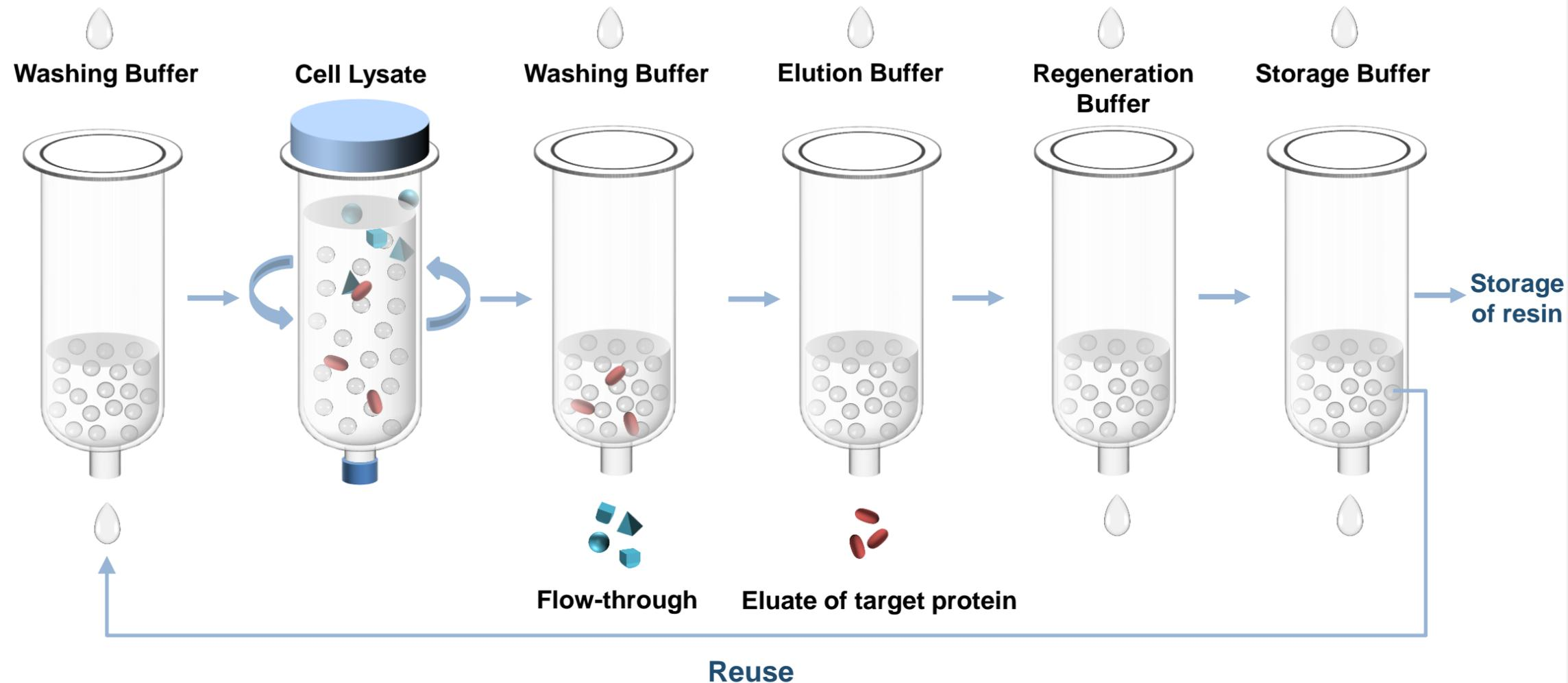
B. Sample Loading: Load sample by incubation or gravity flow.

C. Washing: Wash with 10 bed volumes washing buffer to remove impurities.

D. Elution: Elute protein with 5–15 bed volumes of elution buffer.

E. Regeneration: Flush resin with 15 bed volumes of fresh regeneration buffer (0.1M NaOH).

F. Neutralization & Cleaning: Wash with 2×10 bed volumes storage buffer. Store resin in storage buffer at 2–8 ° C.



3 Result Analysis

Electrophoresis:

Analyze eluted proteins by SDS-PAGE or Western Blot.

Interpretation:

Determine target protein purity and recovery efficiency based on band intensity and position.

