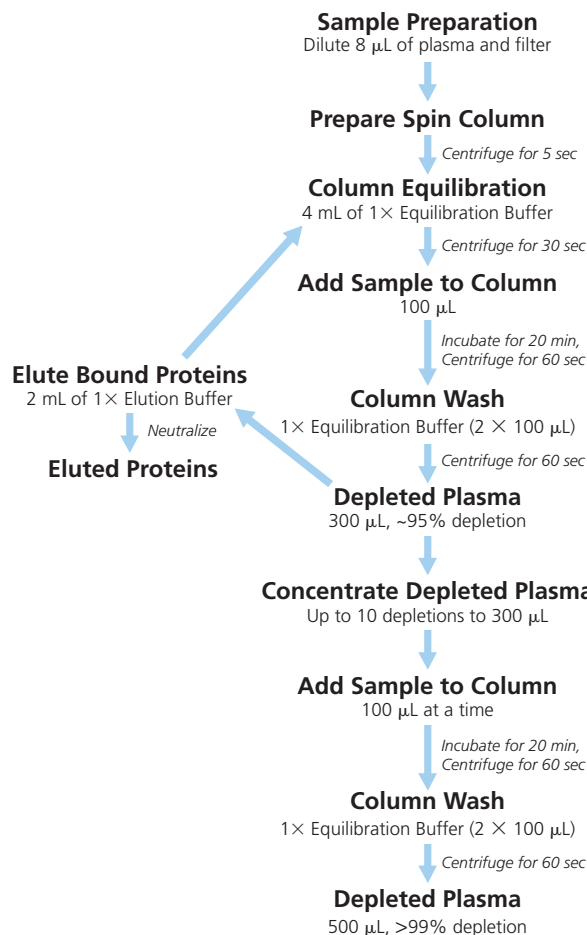


# User Protocol



## Important Reminder

For column storage, the column plug **must** be inserted only after centrifuging the 1 $\times$  Equilibration Buffer from the spin column. Plugging the column without centrifugation could lead to the frit being pushed up by hydraulic pressure.