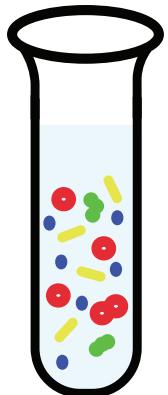
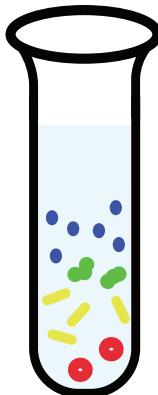


Exosome Isolation: Density Gradient Ultracentrifugation



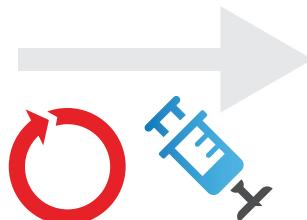
1

Remove cells
and cell debris

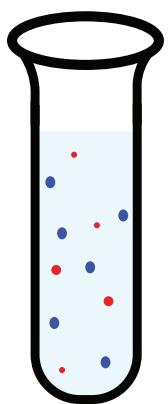


2

Remove smaller debris

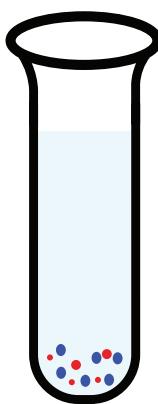


Low force centrifugation:
 $750 \times g$, 15 mins; $2,000 \times g$, 15 mins



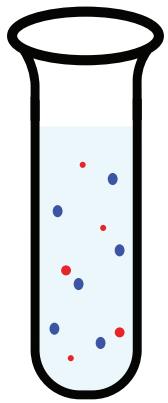
3

Pellet the exosomes



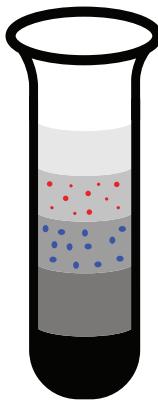
4

Resuspend the pellet



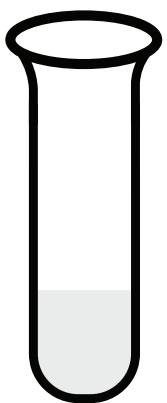
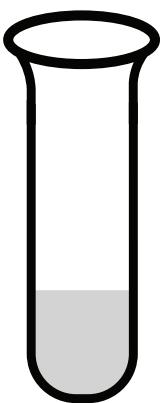
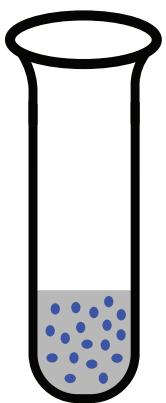
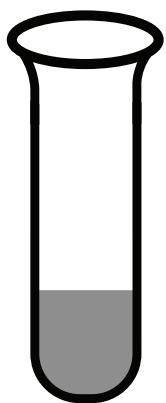
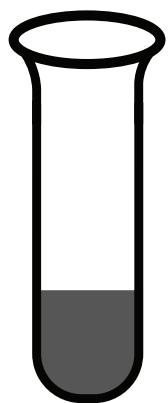
5

Layer onto density gradient
and ultracentrifuge



6

Ultracentrifuge the
gradient fractions



7

Size determination
analysis

8

Experimental
characterization
of exosomes