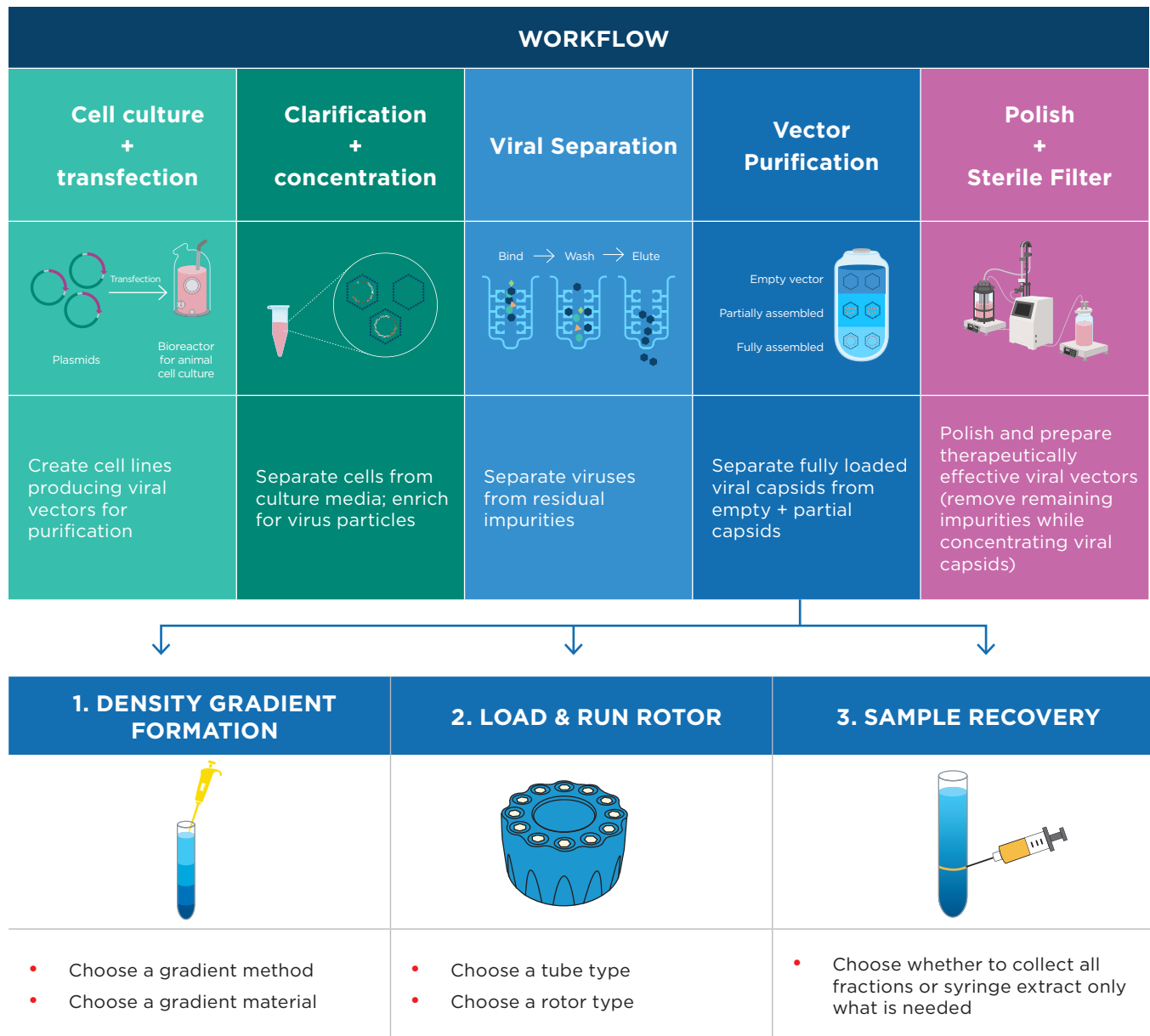




Viral Vector Purification with Ultracentrifugation: How do I separate empty, partial & filled viral capsids?

Vectors such as AAV and adenovirus are powerful delivery tools that are currently used in research, preclinical, and clinical developments. Density gradient ultracentrifugation (DGUC) offers a serotype-independent method of separating empty, intermediate, and full viral particles.

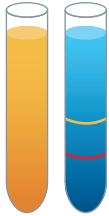
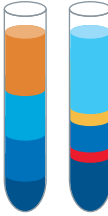
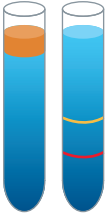




Use Density Gradient Ultracentrifugation to achieve >99% full viral capsids*

* See doi: [10.1089/hgtb.2015.051](https://doi.org/10.1089/hgtb.2015.051) for an example protocol with ultracentrifugation



Considerations in Density Gradient Ultracentrifugation

1. DENSITY GRADIENT FORMATION	GRADIENT METHOD		
	<p>Isopycnic:</p> <ul style="list-style-type: none"> Materials separate based on buoyant density in a self-forming density gradient that results in a continuous density gradient Ideal for separating same size or mass but different buoyant density materials (e.g., empty vs loaded virus) Highest resolution because density gradient is a continuum 	<p>Equilibrium zonal:</p> <ul style="list-style-type: none"> Materials separate by density in a pre-formed density gradient that does not achieve a continuous gradient, but rather has discrete segments of different densities Provides a balance of speed (because the gradient is pre-formed) + purity Less maximum resolving power than isopycnic 	<p>Rate zonal:</p> <ul style="list-style-type: none"> Materials separate based on sedimentation coefficient (e.g., including size and mass) in a pre-formed density gradient Ideal for separating a protein complex from other proteins 
	GRADIENT MATERIAL		
2. LOAD + RUN ROTOR	<p>Iodixanol</p> <ul style="list-style-type: none"> Typically faster protocols Minimal interference with subsequent analysis / analytics Typically used for equilibrium zonal gradients Iodixanol equilibrium zonal density gradients are often selected for rapid, high throughput purification of vectors in discovery or development 	<p>Cesium Chloride (CsCl)</p> <ul style="list-style-type: none"> Better purity and resolution of empty, full, and partial capsids Typically more time-consuming protocol May require lengthier buffer exchange prior to analysis Typically used for isopycnic gradients CsCl isopycnic density gradients are often selected for the highest purity results in manufacturing or when producing reference standards 	
	TUBE SEALING METHOD		
	<p>OptiSeal</p> <ul style="list-style-type: none"> Plug-based seal Faster to seal No physical guarantee of tube remaining sealed during transport / handling 	<p>Quick-Seal</p> <ul style="list-style-type: none"> Permanent, guaranteed seal Heat-based seal Slower to seal 	
TUBE MATERIAL			
<p>Ultra-Clear</p> <ul style="list-style-type: none"> Highly transparent tubes, easier to see bands Not autoclavable, but can be cold sterilized 	<p>Polypropylene</p> <ul style="list-style-type: none"> Resistant to more chemicals than Ultra-Clear tubes Autoclavable Available for Optiseal and Quick-Seal tubes 	<p>Sterile + Certified Free* Tubes</p> <ul style="list-style-type: none"> Available for both Ultra-Clear and polypropylene tubes Only Quick-Seal and open top tube types available 	

*Based on sample results below detectable limits

ROTOR TYPE

2. LOAD + RUN ROTOR (Continue)

Vertical:

- Most efficient rotor for (a) equilibrium zonal separations and (b) isopycnic separations with minimal contaminants that pellet



Near Vertical:

- Ideal for isopycnic separations with crude sample containing significant amounts of contaminant



Fixed Angle:

- Ideal for labs either budget- or space-constrained requiring 1 rotor suitable for both density gradient and pellet separations



Swinging bucket:

- Ideal for rate zonal separations
- Inefficient for other density gradients (e.g., viral separation)

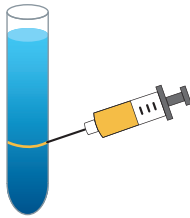


SAMPLE RECOVERY

3. RECOVER SAMPLE

Syringe extraction:

- Precise method to remove only the band of piercing the tube with a syringe to collect the band
- Ideal for well-developed processes where the desired band of interest is known visually or by location.



Fractionation (from middle or bottom of tube):

- Pierce tube with a needle to drain out some of all of the solution in small increments, allowing independent analysis of each fraction.
- Ideal when bands are not visible or when seeking to characterize the composition of multiple bands

