



## OptiMATE Gradient Maker Frequently Asked Questions (FAQs)

### Dispensing FAQs

**1. What should I do if the instrument won't dispense, seems to dispense slowly, or the liquid level doesn't match the dispense volume?**

- Check that the tubing is properly clamped in the pumps.
- Check for the correct tubing connection per the method being used and ensure that the tube is dipped such that the opening is below the liquid level.
- Ensure tubing lines that are in use are not pinched at the manifold.
- Ensure Luer locks are completely connected.

**2. Why do I see bubbles on the top of the tube even after the tube is filled and sealed?**

Bubbles can occur because of the sample composition. Small bubbles are acceptable and should not impact the centrifugation run. However, do not leave a large air space – too much air can cause tube deformation.

**3. Can I leave the tubing in the pumps and for how long?**

Yes, you can leave the tubings in the pumps, typically overnight clamping should not result in any issues. It is strongly recommended to clean the tubing prior to a shutdown of the instrument or longer periods of non-use.

**4. Do I need to disconnect and remove reagent containers after every run?**

- Leaving the dip tube / tubing set in your solutions can cause fouling and reduce the usable life.
- There is no need to disconnect when doing back-to-back runs. However, we recommend disconnecting reagents and sealing with the Luer caps provided and storing appropriately at the end of the day.

**5. How do I avoid bubbles in my tubing?**

- Ensure the dip tube is fully submerged, each reagent bottle has sufficient volume, and all Luer connections are securely attached.
- Ensure tubing set is properly primed and prime again as needed or until bubbles are removed.
- Small, immobile bubbles are generally fine and will not impact your gradient.

**6. Do I need to prime the tubing again if I'm running the same method again?**

No, you don't need to re-prime for back-to-back runs.

**7. How often do I need to calibrate the instrument pumps?**

- Calibrate the pumps at the start of each run.
- If running the same method consecutively, recalibration is not necessary.
- Always recalibrate when switching methods or reagents, or after any tubing has been disconnected.

**8. How often should I replace the probe guide and how can I clean it?**

Clean the probe after every use to avoid chemical deposits. For cleaning, refer to [Document IN-175 Chemical Resistances for Beckman Coulter Centrifugation Products](#).

**9. Can I use third-party tubing?**

No, OptiMATE is only compatible with the tubing provided by Beckman Coulter Life Sciences.

**10. Do I need to weigh tubes before centrifugation if I am loading the sample manually after the OptiMATE instrument dispenses into the tube?**

If all samples loaded are of the exact same volume and gradient profile, then weighing is not required. However, re-calibration is necessary whenever the tubing set is replaced.

**11. How often do I need to change the dispense probe?**

Ensure the dispense probe is thoroughly clean between uses and the connection between the manifold is tight. There is no specification or recommendation for the frequency of probe replacement, as customer workflows and standard operating procedures vary dramatically. Ensure the dispense probe is not clogged or bent.

**12. What density gradient materials can be used with the system?**

Cesium chloride, iodixanol, and sucrose

**13. Are the tubing sets reusable?**

Yes, tubing lines are reusable. They have limited reuse including two cleaning cycles with specified cleaning agents using 1 M sodium hydroxide, and 70% isopropyl alcohol. See instrument IFU for more information.

**14. Can the tubing sets be sterilized?**

Yes, tubing sets can be sterilized via autoclave or gamma irradiation but multiple cycles will reduce the operating lifespan.

## Sealing FAQs

### 15. Can I use third-party tubes?

No, the OptiMATE system is only compatible with ultracentrifuge tubes provided by Beckman Coulter Life Sciences.

### 16. Do I need to remove the seal former when not in use?

A functional seal former does not need to be removed when not in use. However, it should be periodically inspected for residue or deformations and replaced as needed.

### 17. How often should I replace the seal former and how can I clean it?

Ensure the seal former is free of any residue for proper sealing. For cleaning, refer to [Document IN-175 Chemical Resistances for Beckman Coulter Centrifugation Products](#).

### 18. Why didn't the tube seal correctly?

- Ensure the seal former is free of residue and any defects, and replace if needed.
- Ensure the correct tube part number is chosen in the method.

### 19. Why did my tubes from the OptiMATE instrument deform or leak after centrifugation?

- Tubes filled and sealed using the OptiMATE system should perform as expected in the ultracentrifuge. Check the relevant rotor and tube instructions for use (IFU) to ensure the ultracentrifuge run settings were suitable.
- However, some users load their sample manually after gradient dispense, typically if they are loading different samples in a single run. If samples are topped off manually, wick off any excess liquid in the neck of the tube using a laboratory wipe prior to sealing.
- Ensure the seal former is free of residue and any defects, and replace if needed.

### 20. Do I need to squeeze the tubes to check for leaks before centrifugation?

Tubes filled and sealed by the OptiMATE system should not need to be squeezed. Squeezing can disturb the gradient.

## User Interface and Software FAQs

### 21. How do I save a method?

The OptiMATE instrument autosaves as changes are being made. However, published methods cannot be altered. They must be duplicated and renamed before editing.

### 22. How do I add a section to the top of the tube?

Refer to the OptiMATE system IFU. Briefly, add a section to the bottom of the tube, then click

Move Up until it gets to the desired location

### 23. Can I use the same method with different tube types or part numbers?

No, you must create a new method when changing tube types (unless changing only the number of tubes).

#### 24. Why does the tube type matter for a method?

Filling and sealing performance are impacted by tube material, tube size, and tube volume.

#### 25. Can I use the predefined reagents and methods provided with instrument?

Yes, but you should ensure they match your application needs.

#### 26. Do I need to fill tubes all the way to the top?

Check the relevant rotor and tube instructions for use (IFU). Generally, all tubes should be filled to max volume for proper support during centrifugation, except for certain Open-Top, thickwall tubes which can be partially filled.

#### 27. Can I unpublish or edit a published method?

No, once a method is published it is locked and can no longer be edited. It can be duplicated and then edited. It is also possible to delete methods.

For more information and support, please visit our  
OptiMATE instrument and reagents webpages

### OptiMATE

Gradient Maker



<https://becls.co/Optimate>

### OptiMATE

Density Gradient Reagents



<https://becls.co/OptimateReagents>