

Ksep® 50



Ksep® 50 Product Tutorial



<https://sar.to/ksep50>



https://youtu.be/yUKAE7_CUHI



In this video series, we'll take you through the installation of the chamber and tube sets for the Ksep® 50 centrifugation system, and show you how to calibrate the pump, test the bubble sensors, start the system, create a recipe and export data.

