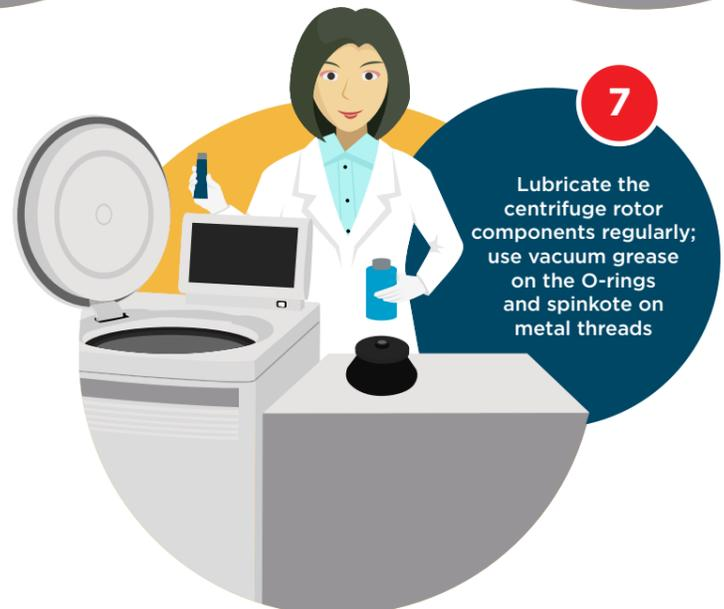
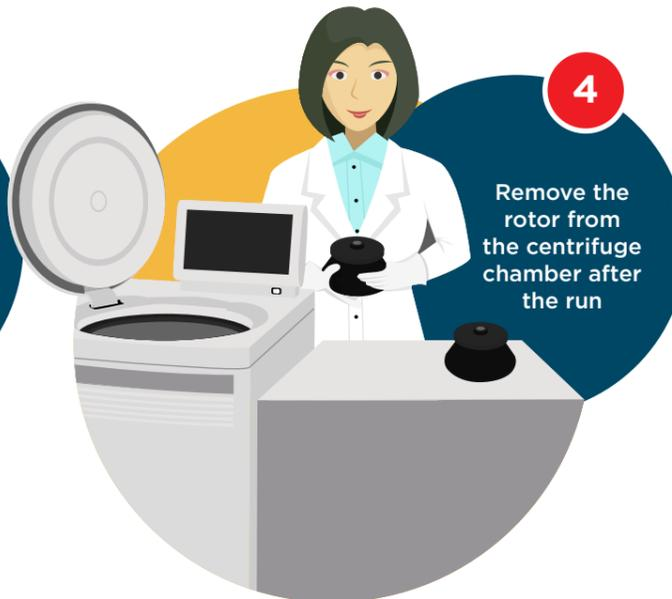
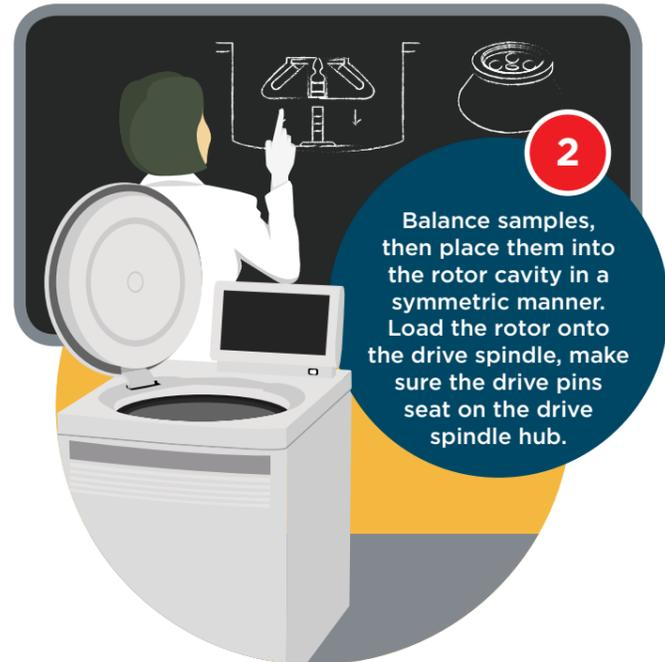




Take good care of your instrument Start with the basics

Operating tips for Beckman Coulter Avanti High Performance Centrifuges



Beckman Coulter
Life Sciences

 <http://www.beckman.com/contact-us>



© 2016 Beckman Coulter, Inc. All rights reserved. Beckman Coulter, the stylized logo, and the Beckman Coulter product and service marks mentioned herein are trademarks or registered trademarks of Beckman Coulter, Inc. in the United States and other countries.

For Beckman Coulter's worldwide office locations and phone numbers, please visit "Contact Us" at beckman.com