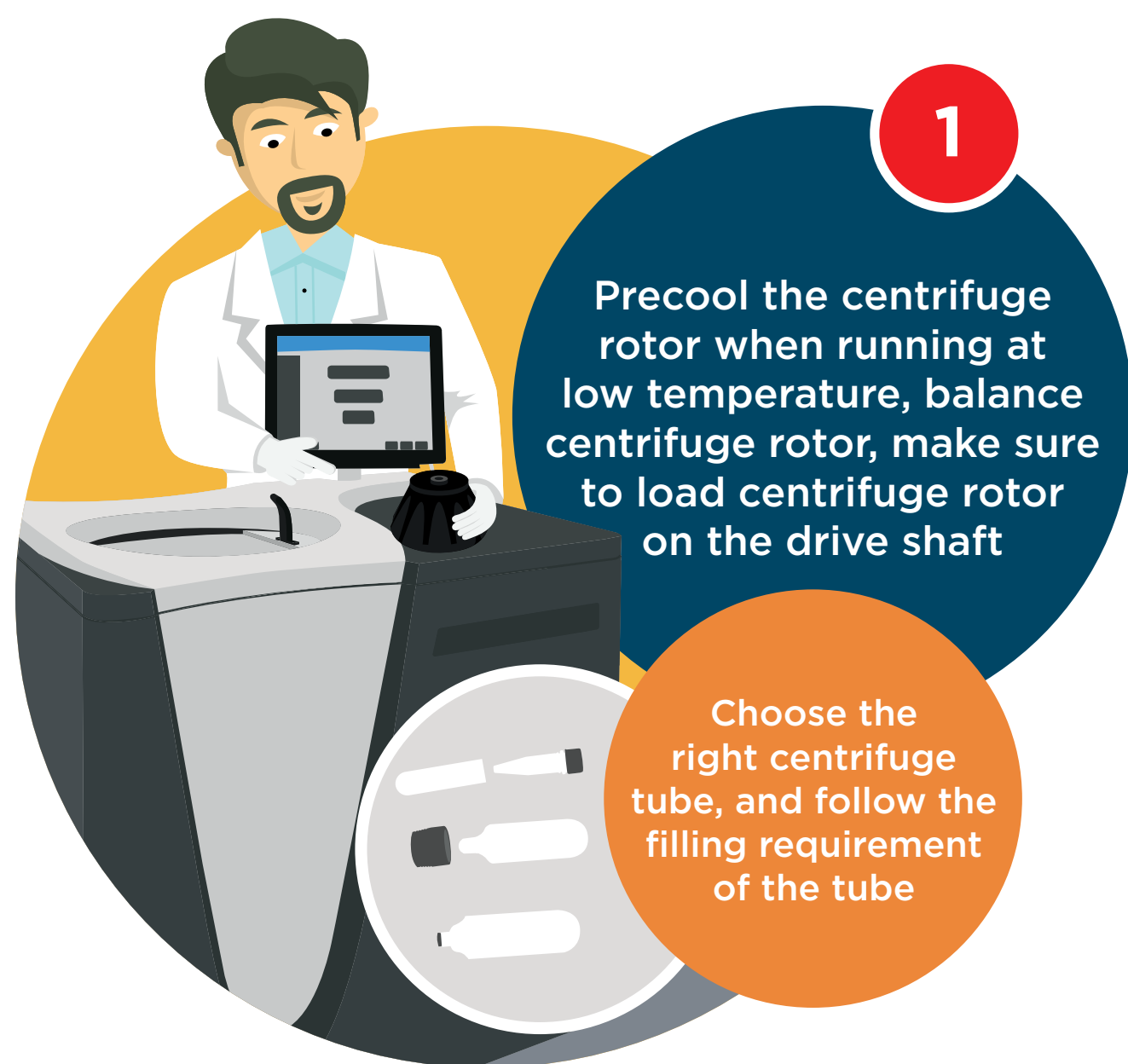




# Take good care of your instrument Start with the basics

Operating tips for a Beckman Coulter Ultracentrifuge



**1**


Precool the centrifuge rotor when running at low temperature, balance centrifuge rotor, make sure to load centrifuge rotor on the drive shaft

Choose the right centrifuge tube, and follow the filling requirement of the tube



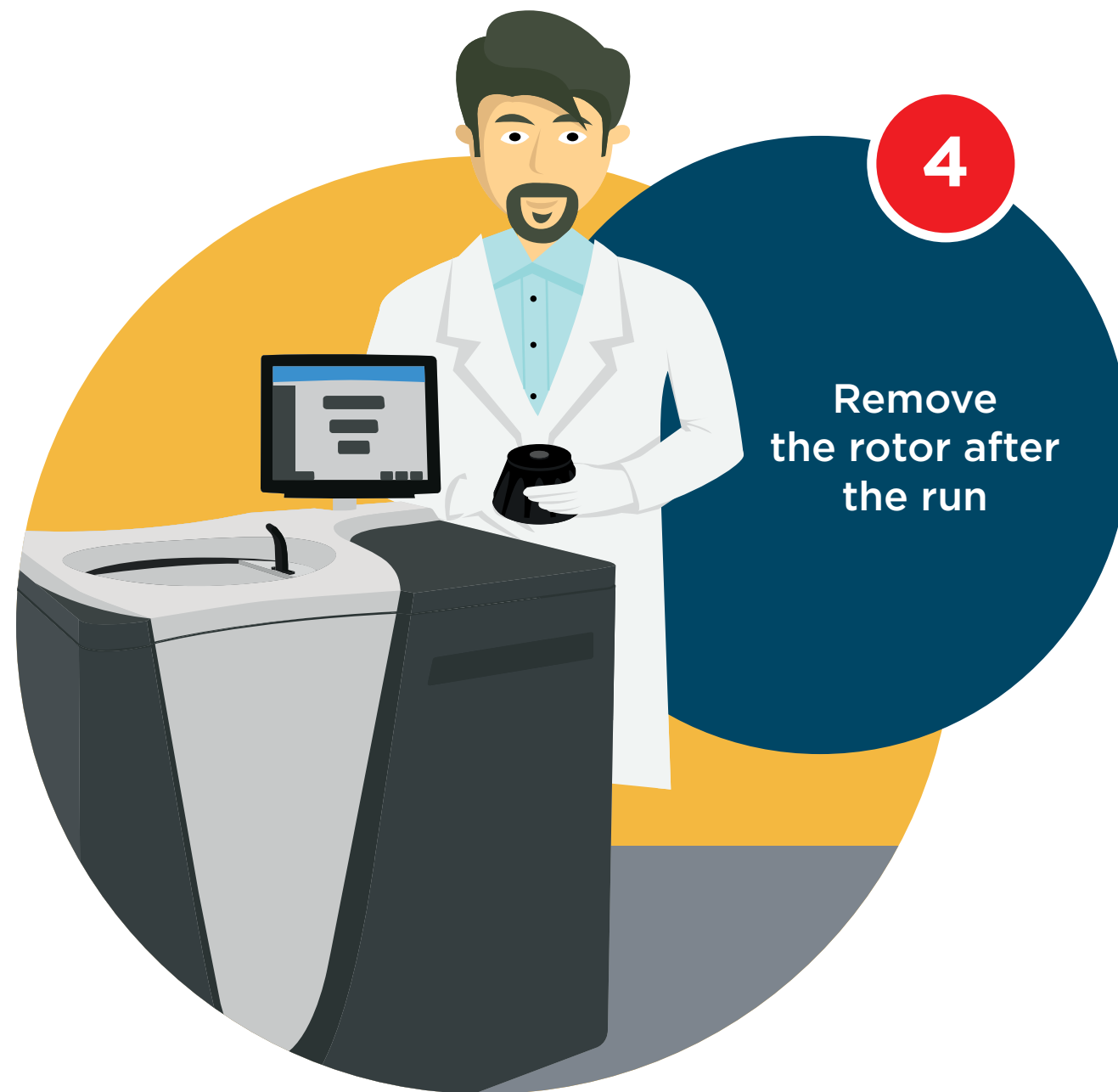
**2**

When using the swinging bucket rotor, match numbered caps with numbered buckets and screw the caps in until there is metal-to-metal contact



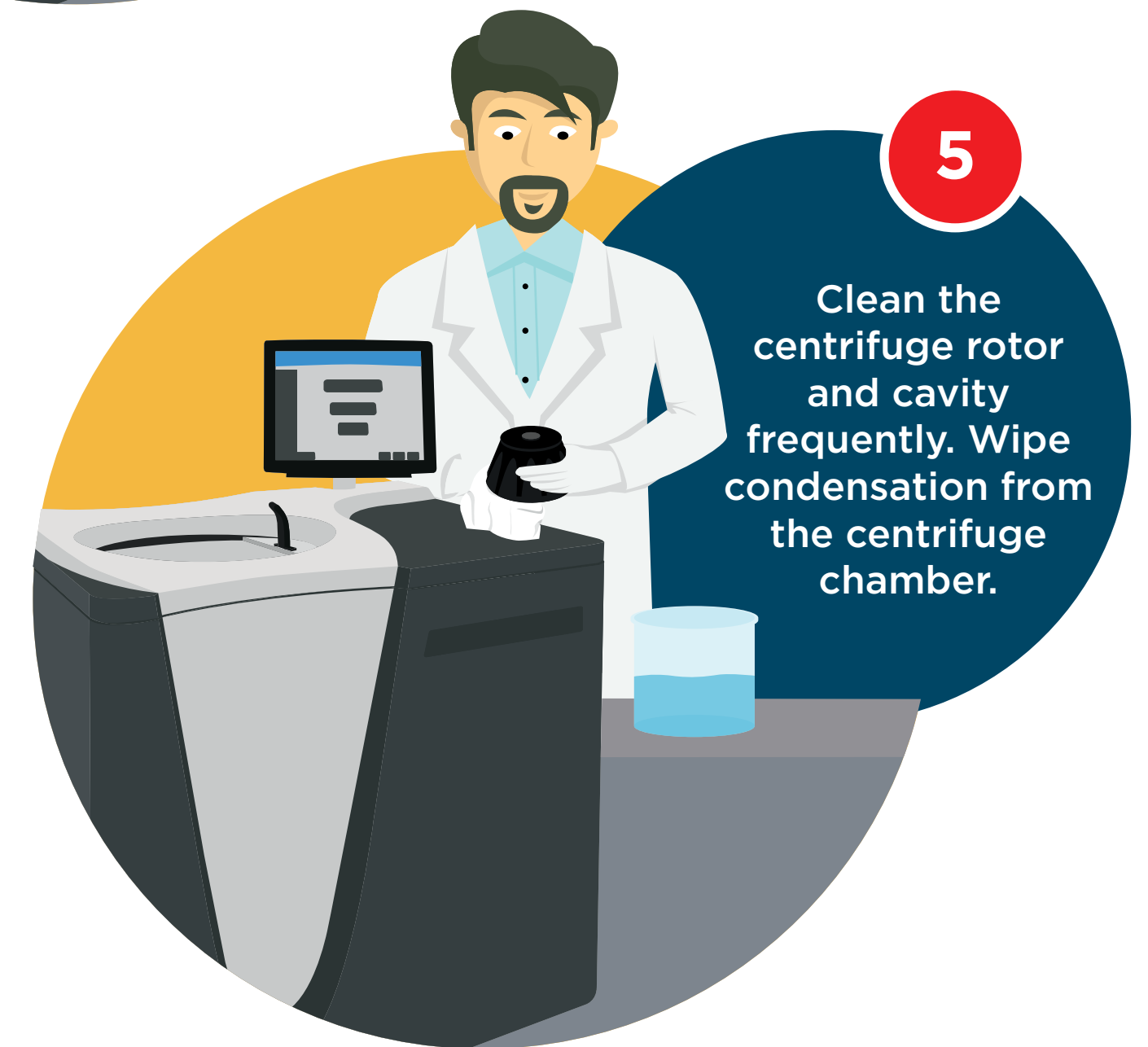
**3**

Never use a sharp object on the touch screen panel to set parameters



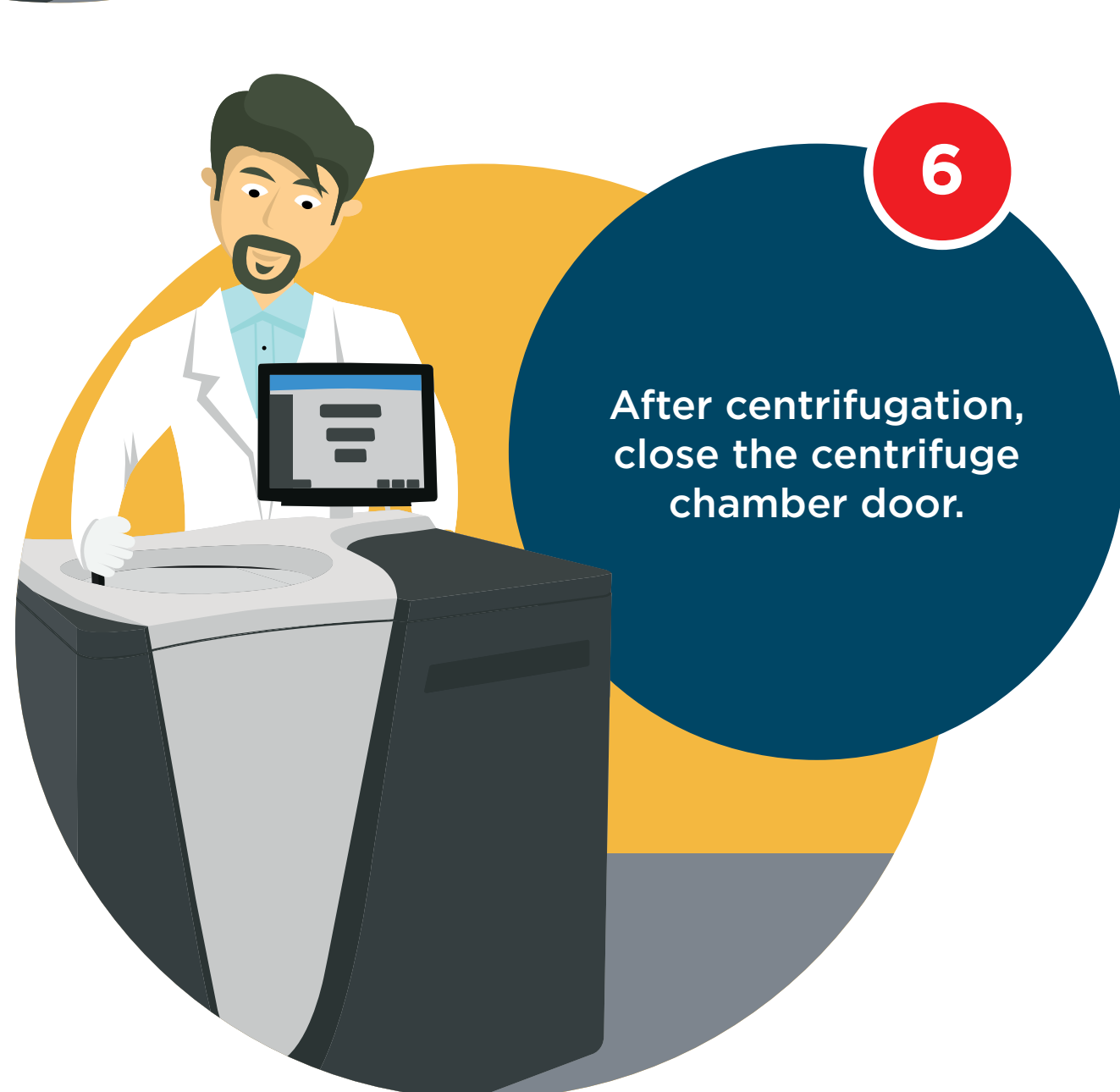
**4**

Remove the rotor after the run



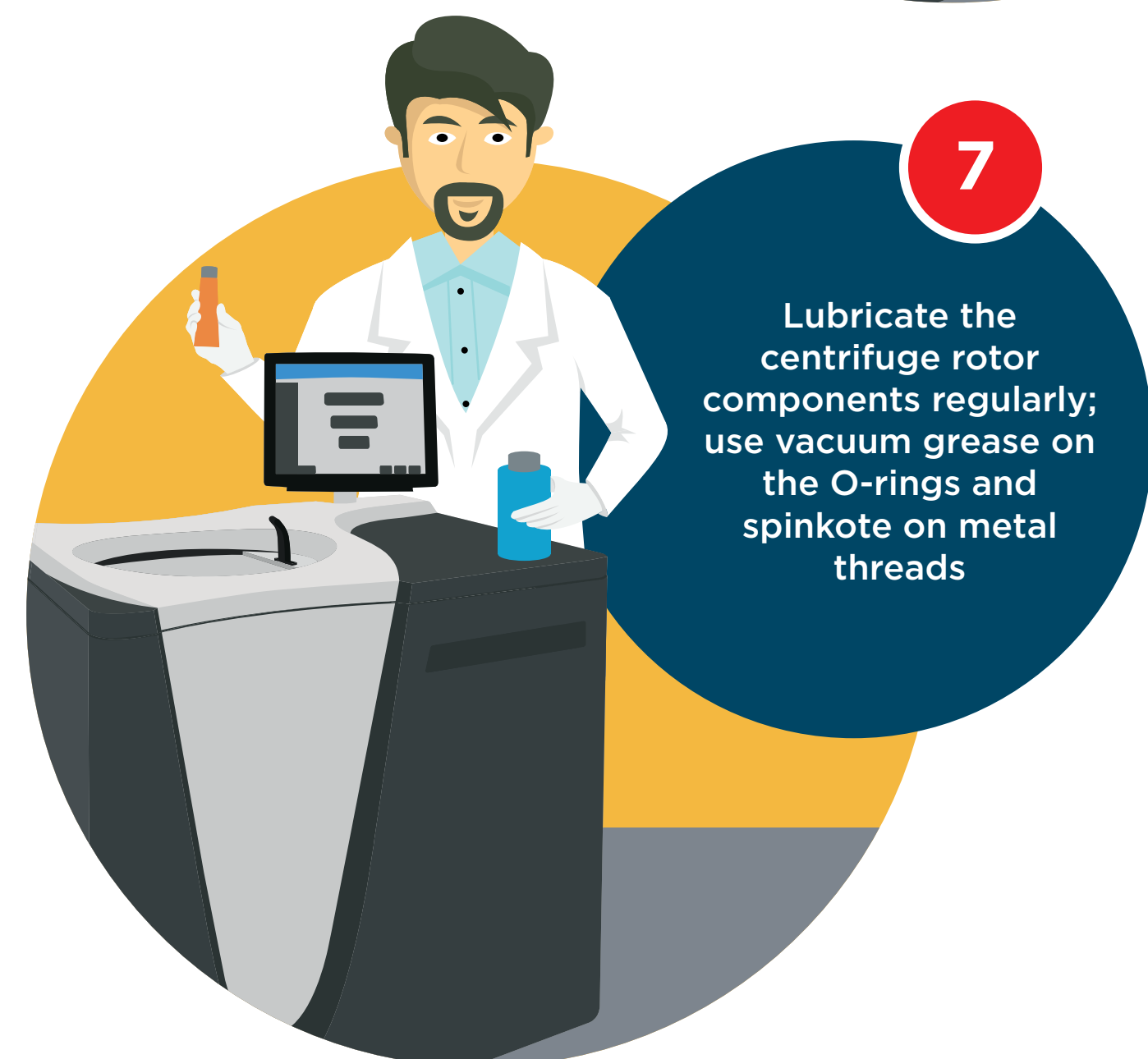
**5**

Clean the centrifuge rotor and cavity frequently. Wipe condensation from the centrifuge chamber.



**6**

After centrifugation, close the centrifuge chamber door.



**7**

Lubricate the centrifuge rotor components regularly; use vacuum grease on the O-rings and spinkote on metal threads

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](http://beckman.com)