

Microplate Centrifuge Unpacking Guide

Before you start

Move the shipping box to the area where you will be installing the device. Make sure two people are available to unpack the device.

WARNING The fully packed Microplate Centrifuge weighs approximately 33 kg (72.8 lb). The Centrifuge device weighs 30 kg (66 lb). Attempting to move the device without assistance could cause personal injury. Request assistance and use proper lifting techniques when lifting the device.

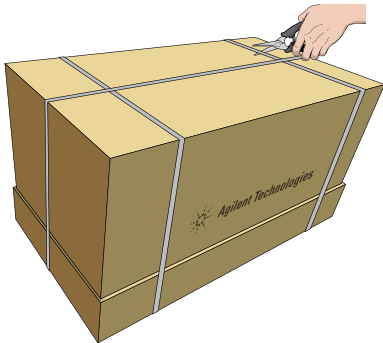
Unpacking the shipping box

IMPORTANT Save the packaging materials in case you need to move or ship the device.

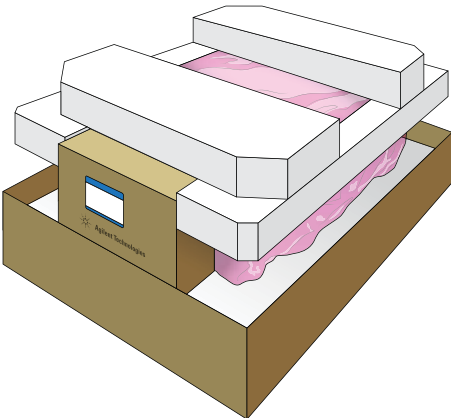
To unpack the Microplate Centrifuge:

- 1 Cut the straps that secure the shipping box.

WARNING The strap will spring back with force when cut. To avoid potential injury, stay clear of the sides of the shipping box when cutting the packing strap.

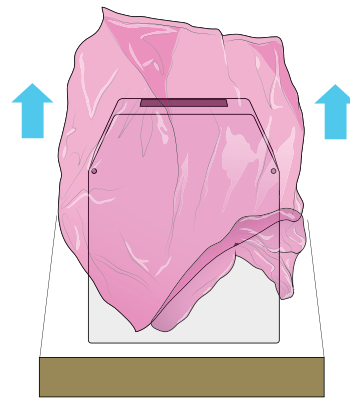


- 2 Lift the top cardboard cover off of the box, and then remove the top foam packing insert.



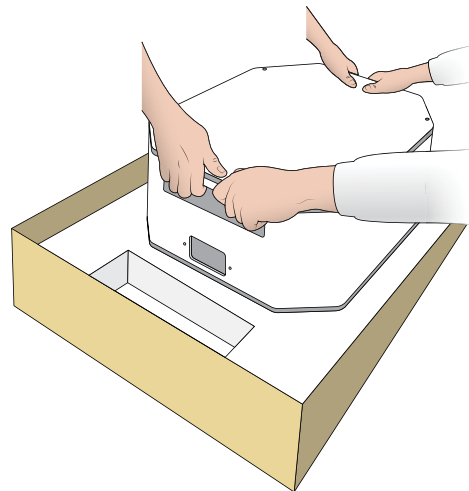
- 3 Remove the small box, and set it aside temporarily.
- 4 Remove the top foam packing insert from the device.
- 5 While the Centrifuge is still in the box, remove the Important Documents envelope from the plastic bag. Set aside the envelope temporarily.
- 6 While the Centrifuge is still resting on the packing foam in the shipping box, use the door opening to carefully tilt the device up vertically. The rear end should remain resting on the foam and the closed end of the bag should be facing up.

While supporting the Centrifuge to prevent it from falling, lift the plastic bag off of the device.



- 7 Gently, lower the Centrifuge down onto the foam again.
- 8 Refer to the following figure for the correct lift points on the Centrifuge.

Figure Correct lift points on the Centrifuge



9 With a person positioned on each side of the device, grasp the Microplate Centrifuge using the front door opening and the rear connection panel recess as lift points.

While firmly holding the device, carefully lift it out of the shipping container and onto the laboratory table.

10 Open the small box that you set aside in [step 3](#). Carefully, unwrap the packing foam and remove the pair of centrifuge buckets.

Inspecting the contents

If an item is defective on arrival, contact Agilent Technologies as soon as possible.

- In North America, call 1.800.227.9770.
- For other locations, contact your region's Agilent Technologies office:
<http://www.chem.agilent.com/en-US/ContactUS/Pages/ContactUs.aspx>.

Note: The power and communication cables, air fittings, and other items are shipped separately.

About installation, setup, and operation

For safety information, laboratory setup requirements, installation instructions, and operating instructions see the *Microplate Centrifuge User Guide*.

For operating instructions, see the *Microplate Centrifuge Quick Guide*.

The user guides are available from the Help menu of the VWorks software or as PDF files on the software CD. For the latest versions, go to www.agilent.com/chem/askb.

Contacting Agilent Technologies

Technical Support: 1.800.979.4811 (US only) or +1.408.345.8011 (rest of the world)
service.automation@agilent.com

Customer Service: 1.866.428.9811 or +1.408.345.8356
customerservice.automation@agilent.com

European Service: +44 (0)8457125292
euroservice.automation@agilent.com

Documentation feedback:
documentation.automation@agilent.com

Web: www.agilent.com/lifesciences/automation