

BenchBot Robot Unpacking Guide

Before you start

The BenchBot Robot is shipped in the following boxes:

- *BenchBot Robot box.* Contains the robot only.
- *Accessories box.* Contains the emergency-stop pendant (standalone/component configurations only), cables, teaching plate, regrip station, screws, software CD, and user documentation.

Note: The integration plates and the safety equipment supplied with the workstation configuration are shipped in additional boxes. The size and number of these boxes depend on the size of the workstation.

WARNING The robot shipping box, containing the BenchBot Robot, with pallet weighs 35 kg (77 lb). Use a small pallet jack to lift and move the box.

Move the shipping boxes to the area where you will be installing the robot.

WARNING The BenchBot Robot weighs 19 kg (42 lb). Do not attempt to move the robot without assistance. Use proper techniques when lifting the robot.

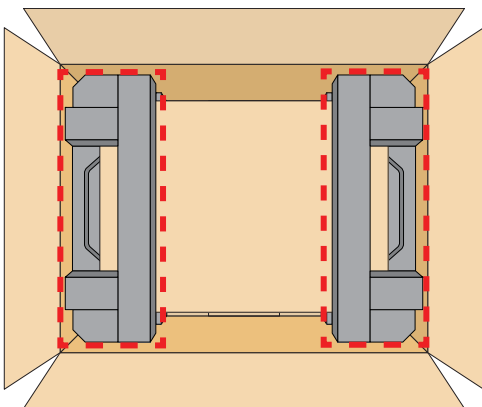
Make sure two people are available to lift and move the robot from the box.

Unpacking the shipping boxes

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

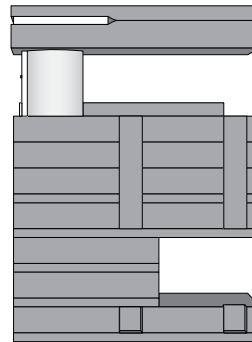
To unpack the BenchBot Robot box:

- 1 Cut and remove the plastic straps that hold the shipping box to the pallet.
- 2 Open the shipping box, and remove the two foam blocks covering the smaller box (box 2) inside.

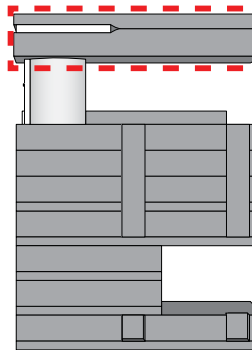


- 3 With two people, remove box 2, place it on the floor, and lift open the lid.
- 4 With two people, carefully lift the robot out of box 2 and place the robot upside-down on a flat, stable surface where it will be possible for you to access all sides of the robot.

Figure Robot wrapped in foam blocks, side view, upside down

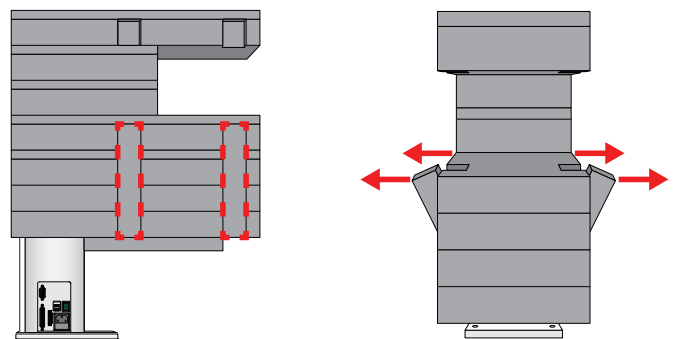


- 5 Remove the two foam blocks from the base.

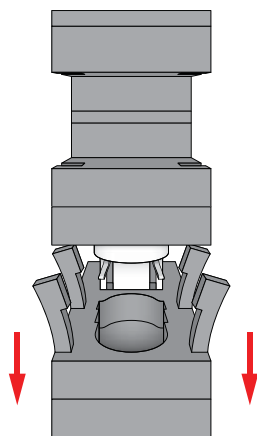


- 6 Turn the robot right-side up so that it is sitting on its base with the front side facing you, and then pull open the top of the four foam tabs as shown.

Figure Robot on its base, side view (left) and front view (right)



- 7 Remove the large foam block by pulling it downward and toward the front of the robot.



- 8 Remove the top foam block.
- 9 Unpack the accessories box and, if applicable, the boxes containing the integration plates and the safety equipment.

Inspecting the contents

Inspect all items for completeness and potential shipping damage. If an item is defective on arrival, contact the Automation Solutions Business Center 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>.

After you unpack the BenchBot Robot box and the accessories box, make sure you have the following:

Part name	Part number
BenchBot Robot	G5486-00001
M6 x 16 screws (4), robot mounting	G5550-02702
Teaching plate	G5550-23357
Regrip station assembly:	
Regrip station	G5550-20020
FHCS M6 x 35 (1), station mounting	G5550-02696
Gripper pad replacement kit	G5486-60064
Power cable	Request by country
Emergency-stop pendant (accessories box for standalone/component configurations or safety equipment box for workstation configurations)	G5486-60026
Ethernet cable	G5550-09002

Part name	Part number
VWorks software CD:	
Benchtop license, or	08330-402
System license	08330-403
<i>BenchBot Robot Safety and Installation Guide</i>	G5486-90001

About installation, setup, and operation

For safety information, laboratory setup requirements, and installation instructions, see the *BenchBot Robot Safety and Installation Guide*.

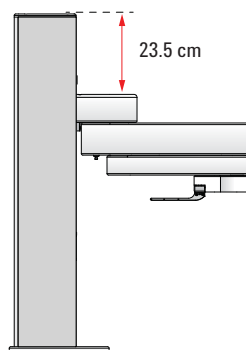
For software setup and operating instructions, see the *BenchBot Robot User Guide*.

Packing for storage and transport

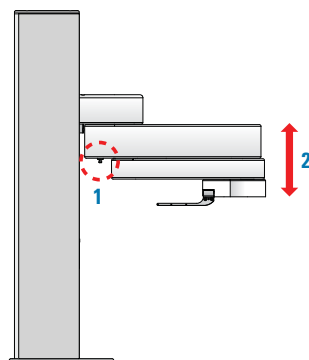
If you need to remove the robot from the attachment surface and package the robot for storage or transport, follow the instructions in this section.

To prepare for packing the robot:

- 1 Home the robot. For instructions, see the *BenchBot Robot User Guide*.
- 2 Check that the top of the shoulder is 23.5 cm (9.25 in) from the top of the mast. This is the correct packing position.



To change the robot position, with the robot power turned on, support the weight of the robot arm and press the z-axis brake release button (1). Raise or lower the arm so that the top of the shoulder is 23.5 cm (9.25 in) from the top of the mast (2).

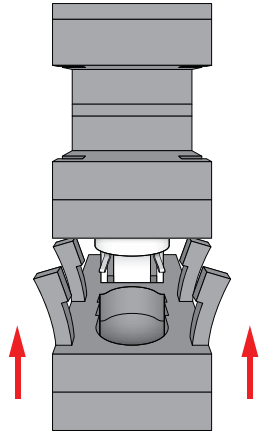


- 3 Turn off the robot and disconnect it from the power source.

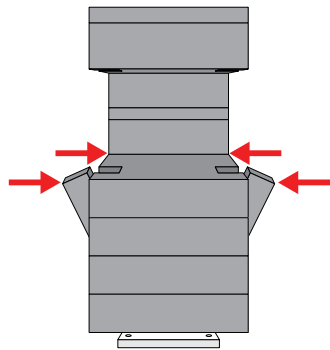
To pack the BenchBot Robot:

- 1 Place the top foam block over the top of the mast.
- 2 Open the four tabs in the large foam block, place the foam block at the base of the robot, and then carefully move it upward so that the robot grippers are positioned within the cutout.

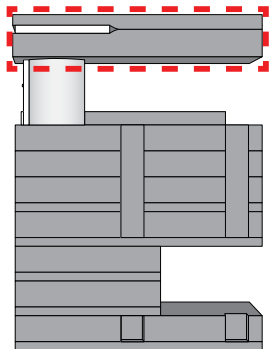
Figure



- 3 Push closed the four tabs in the large foam block.



- 4 Carefully turn the robot upside down.
- 5 Place the thin foam blocks around the base.



- 6 Lift the robot into the smaller box (box 2).
- 7 Place the lid on box 2.
- 8 Place box 2 in the shipping box and place the shipping foam blocks on top of box 2. Close the shipping box and tape it closed.
- 9 Place the shipping box on the pallet, and use plastic straps to secure it.

Be sure to pack the accessories box. Include the teaching plate, regrip station assembly, cables, screws, software CD, and user documentation in the accessories box. Tape the box closed.

If applicable, pack the following:

- *Integration plate boxes.* Include the integration plates and screws.
- *Safety equipment boxes.* Include the BenchBot Disable Hub and all cables, the emergency-stop pendant, the Light Curtain or the shield panels, and all screws and tools required to install the safety equipment.

Accessing product user documentation

Agilent Technologies product user documentation is available in the following locations:

- *Online help.* The product knowledge base is available from the Help menu within the automation software.
- *PDF files.* Product manuals are provided as PDF files on the automation software CD.
- *Agilent Technologies website.* You can search the product knowledge base or download the latest version of any PDF file from the Agilent Technologies website at www.agilent.com/lifesciences/automation.

Contacting Agilent Technologies

Business Center: 1.800.227.9770
(for missing or damaged components)

Technical Support: 1.800.979.4811 or +1.408.345.8011
service.automation@agilent.com

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

European Service: +44 (0)1763850230
euroservice.automation@agilent.com

Documentation feedback:
documentation.automation@agilent.com

Web: www.agilent.com/lifesciences/automation

