

Bravo Platform Unpacking Guide

This guide describes how to unpack the Bravo device from the shipping carton. Depending on the order, the shipment can include additional packages for the Bravo utilities, pipetting head, computer, and accessories.

Before you start

- 1 **Site specifications.** Ensure the installation site meets the requirements described in the [Bravo Automated Liquid-Handling Platform Safety and Installation Guide](#).
- 2 **Tools and equipment.** Ensure that you have a cross-tip screwdriver, 2.5-mm and 3-mm hex wrenches, and a cart for moving the Bravo device.
- 3 **Personnel requirements.** Make sure two people are available to unpack the Bravo device.

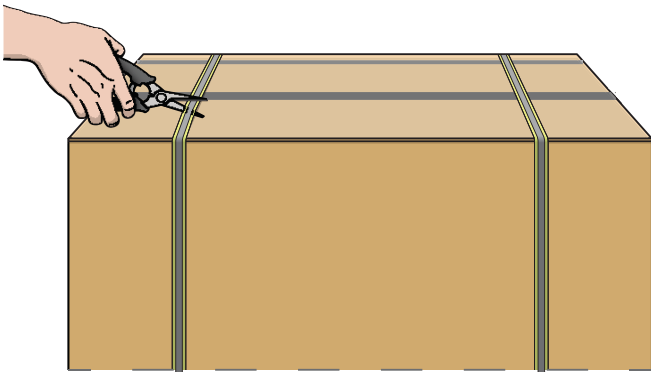
WARNING The Bravo device weighs approximately 52.1 kg (114.9 lb). Attempting to move the device without assistance could cause personal injury. Request assistance and use proper lifting techniques when lifting the Bravo device.

CAUTION Save the packing materials and shipping carton for potential future moves. Packing the Bravo device using other materials might damage the device and void your warranty.

Unpacking the carton

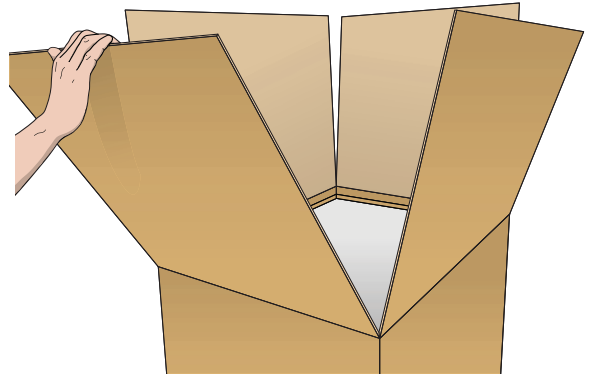
To unpack the Bravo device:

- 1 Cut the straps that secure the Bravo carton.



WARNING Ensure that you stay clear of the sides of the Bravo carton when cutting the packing strap. The strap will spring back with force when cut.

- 2 Open the top flaps of the carton.



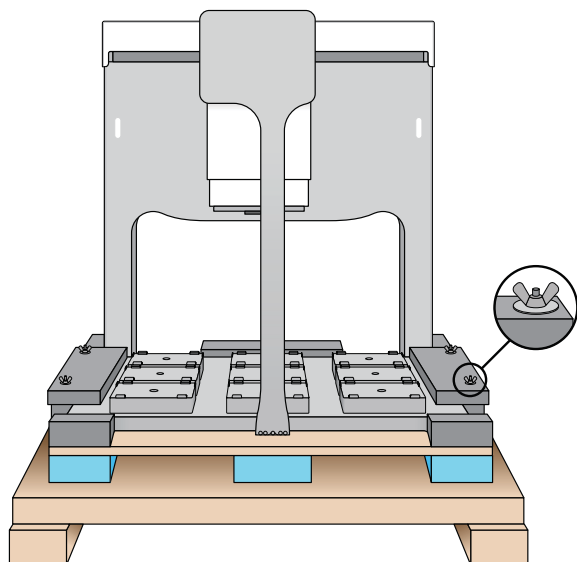
- 3 With two people, one on either side of the carton, grasp the carton at the edges and carefully lift the carton up and off of the Bravo device. The Bravo device remains on the crate base.



Note: The carton interior contains the foam that protects the device during shipping.

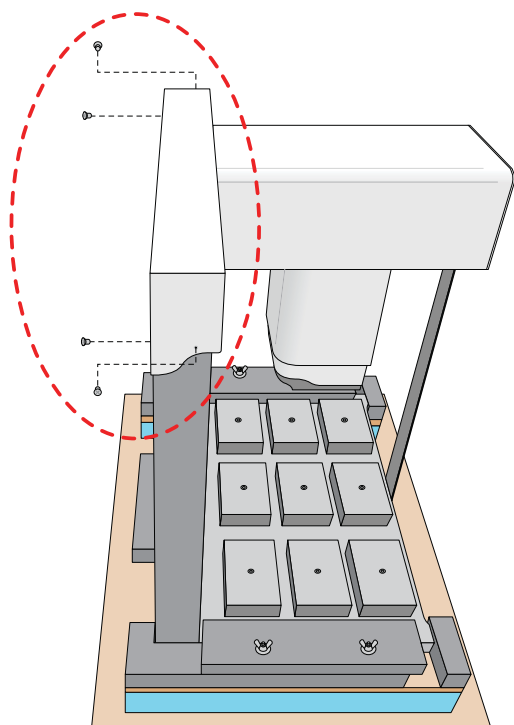


- 4 Remove the plastic covering from the Bravo device.
- 5 Unscrew the four wing nuts and remove the black bars that hold the Bravo deck to the crate base.

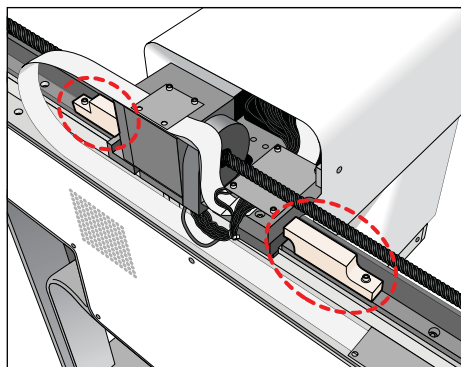


WARNING Do not remove the exterior covers of the Bravo device, except to remove the shipping stops as described in the following steps. Using the device with the exterior covers removed can cause injuries and damage the device.

- 6 Remove the four crosshead screws on the sides and back of the Bravo top cover. Remove the cover.

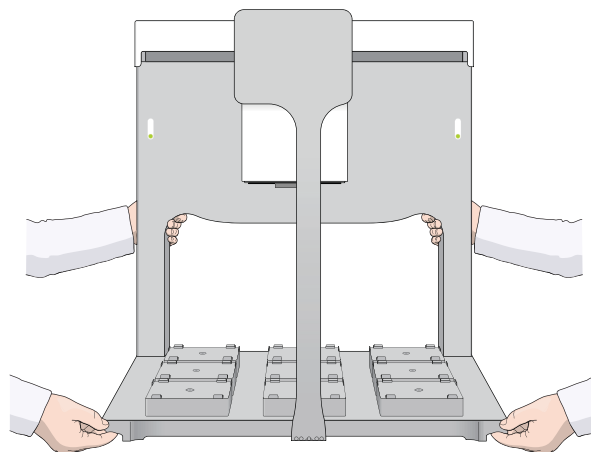


- 7 Using a 3-mm hex wrench, remove the two shipping stops under the top cover.



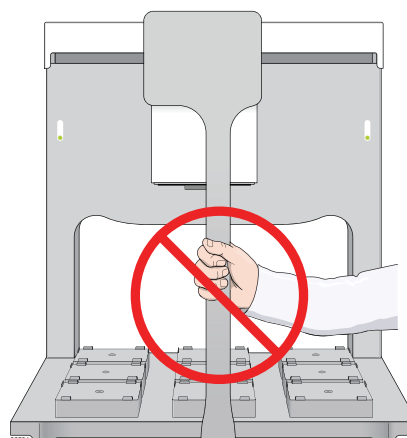
Note: These shipping stops prevent the head mount from moving in the x -axis during shipping.

- 8 Re-attach the Bravo top cover.
- 9 Using the following lift points, lift the Bravo device onto the lab bench or table.

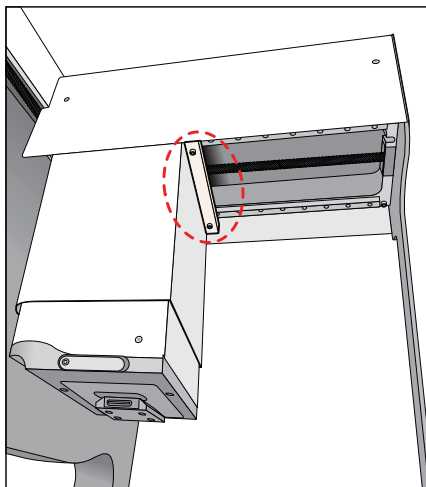


WARNING The Bravo device weighs approximately 52.1 kg (114.9 lb). Attempting to move the device without assistance could cause personal injury. Request assistance and use proper lifting techniques when lifting the Bravo device.

CAUTION Never lift from, or tug on, the tie bar.

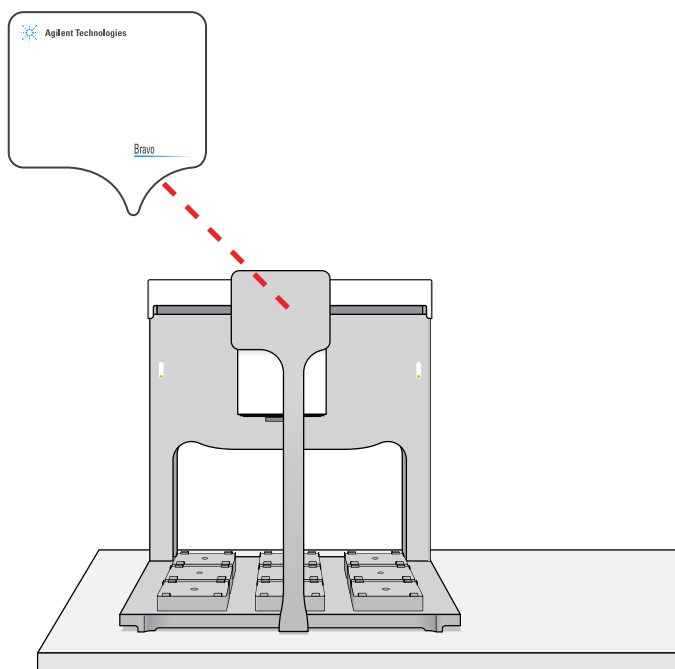


- 10 Using a 2.5-mm hex wrench, remove the shipping stop at the front of the head mount.



Note: This shipping stop prevents the head mount from moving in the y-axis during shipping.

- 11 Position the white face plate, which displays the Agilent logo, at the top of the Bravo tie bar. The face plate is held in place magnetically.



Inspecting the shipped items

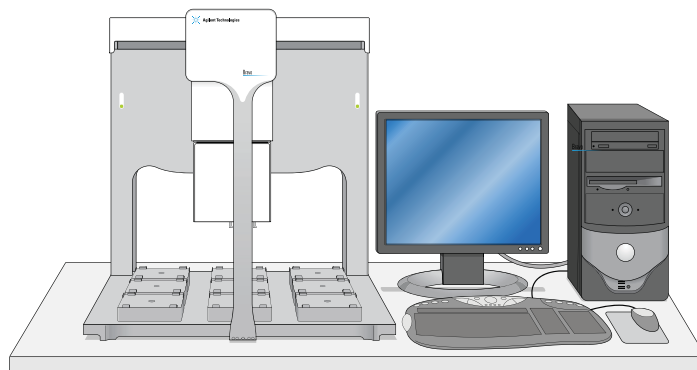
Inspect all items for completeness and shipping damage. If an item is defective on arrival, contact the Automation Solutions Business Center as soon as possible. In North America, call 1.800.227.9770. In other locations, contact your region's Agilent Technologies office: <http://www.chem.agilent.com/en-US/ContactUS/Pages/ContactUs.aspx>

After you unpack the Bravo shipping packages, make sure you have the following items:

- Bravo device
- Pendant, if applicable
- Power, serial, and crossover Ethernet cables
- Pipetting head, if applicable
- Light Curtain and safety shields, if applicable
- Computer, if applicable
- Software CD-ROM
- *Bravo Automated Liquid-Handling Platform Safety and Installation Guide*

About installation

See the *Bravo Automated Liquid-Handling Platform Safety and Installation Guide* for instructions.



Accessing product user documentation

You can access the product Knowledge Base from the help menu within the automation software or online at www.agilent.com/lifesciences/automation.

Contacting Agilent Technologies

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
orders.automation@agilent.com

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

Documentation feedback:
documentation.automation@agilent.com

Web: www.agilent.com/lifesciences/automation