

Automated Centrifuge Loader Unpacking Guide

Before you start

Move the shipping box to the area where you will be installing the Microplate Loader. The fully packed Loader weighs approximately 13.3 kg (29.3 lb).

Ensure that you have the following tools:

- Scissors or other tool for cutting package straps
- 11-mm and 7-mm open wrenches
- Flathead screwdriver
- 3-mm hex wrench

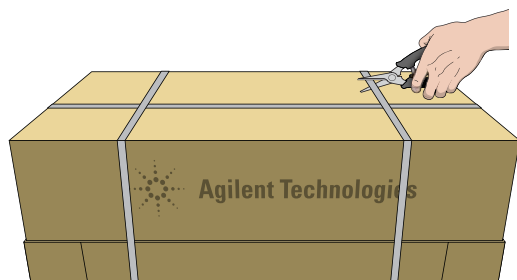
Unpacking the shipping box

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

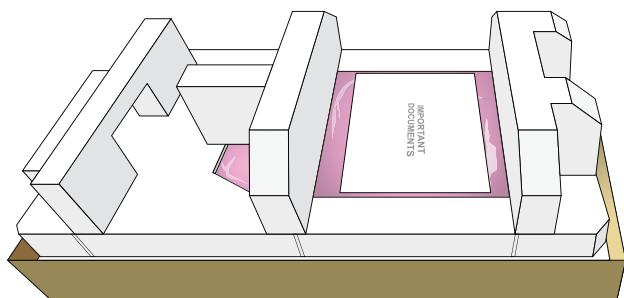
To unpack the Microplate Loader:

- 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.



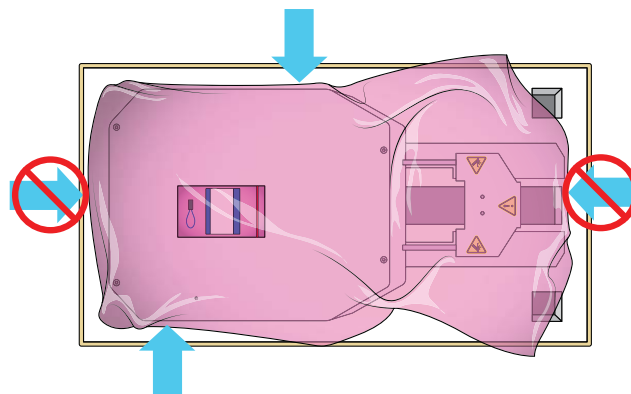
- 2 Lift the top cardboard cover off of the box.
- 3 Remove the Important Documents envelope, which contains the Declaration of Incorporation. Set it aside temporarily.



- 4 Remove the top foam packing insert from the device.
- 5 Ensure that you understand how to lift the Loader using the correct lift points, as the following figure shows.

CAUTION Do not use the front end of the Loader as a lift point. Lifting the Loader by the front end can damage the baseplate, which is easily bent.

Figure Correct and incorrect lift points



- 6 While the device is still resting on the packing foam, carefully, remove the plastic bag that covers the device.

Temporarily, set aside the small bag that contains the e-stop jumper.

Note: Depending on the integration, the jumper might be required instead of an emergency-stop pendant to close the safety interlock circuit.

- 7 With your hands positioned on each side of the device at the correct lift points shown in the preceding figure, reach down between the foam and the device to grasp the baseplate.

While firmly holding the baseplate, lift the Loader onto a laboratory table so that the front end with the Loader shipping bracket is off of the table surface.

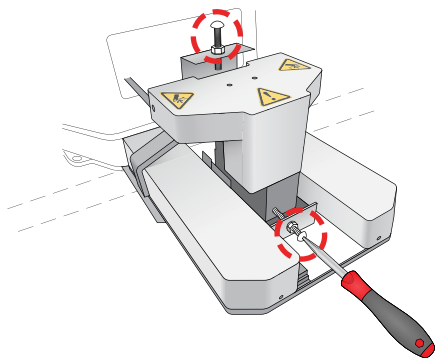
CAUTION Placing the entire device on a flat surface while the screws and the shipping bracket are attached to the Loader will cause potential damage to the device. Ensure that you keep the front Loader shipping bracket suspended off of the table surface until after you uninstall the shipping bracket.



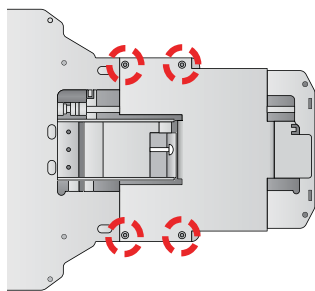
- 8 Uninstall the shipping bracket from the Loader, as follows:

CAUTION Use care to avoid scratching the surfaces of the Loader when uninstalling the shipping bracket and cover.

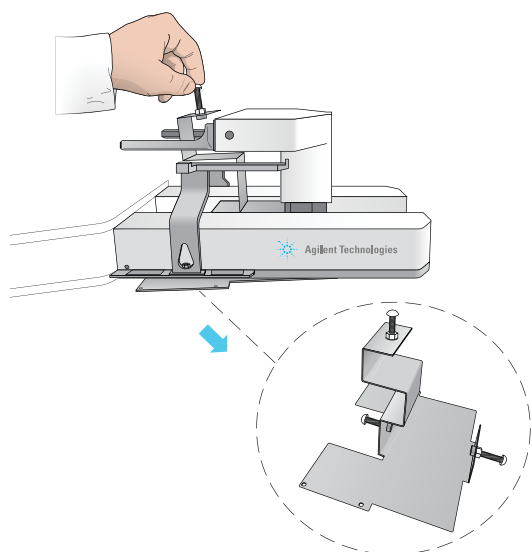
- a To loosen the two nylon fasteners that secure the bracket at the top and front of the Loader, use an 11-mm open wrench to loosen the two nylon nuts, and then use a flathead screwdriver to loosen the nylon screws.



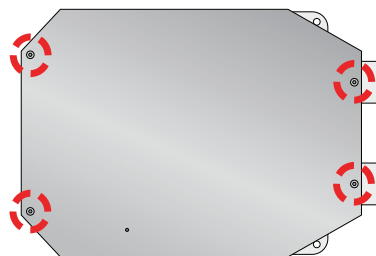
- b Using a 3-mm hex wrench, remove the four socket-head screws that attach the shipping bracket to the bottom surface of the Loader.



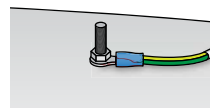
- c Carefully, remove the shipping bracket.



- 9 Remove the four screws that secure the shipping cover to the integration end of the Loader, as the following figure shows.



- 10 Use a 7-mm open wrench to disconnect the ground cable from the shipping cover.



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>.

About installation, setup, and operation

For instructions on how to integrate the Loader with the Centrifuge, contact Automation Solutions Technical Support.

For safety information, laboratory setup requirements, and installation instructions for the Centrifuge with Loader, see the *Automated Centrifuge Loader User Guide*.

For operating instructions, see the *Centrifuge with Loader 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 VWorks Knowledge Base at 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