

Microplate Exchanger Unpacking Guide

Before you start

The Microplate Exchanger is shipped in three cardboard boxes. The largest box contains the Microplate Exchanger, the medium-sized box contains the power supply, and the small box contains the power cord.

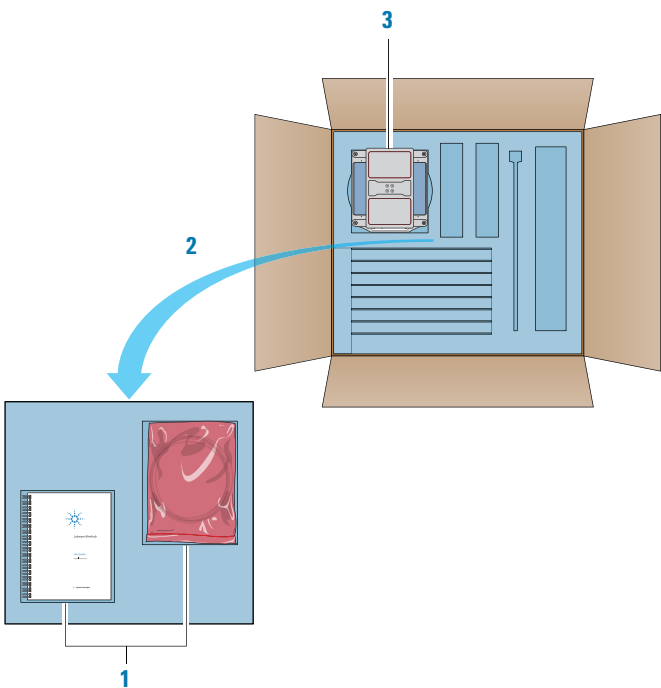
WARNING The Microplate Exchanger shipping box weighs 18 kg (39 lb). Make sure at least two people are available to lift and move the box. Attempting to move the box without assistance could cause personal injury. Use proper techniques when lifting the box.

Move the shipping boxes to the area where you will be installing the device. Leave the boxes on the ground to facilitate the unpacking process.

Unpacking the shipping box

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

To unpack the Microplate Exchanger:



- 1 Open the largest box, and then remove the user documentation and the plastic bag containing cables and screws from the foam block.
- 2 Remove the foam block.

- 3 Carefully lift the Microplate Exchanger base out of the box.
- 4 Open the medium-sized shipping box and remove the power supply from the box.
- 5 Open the small shipping box and remove the power cord from the box.

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 shipping boxes, make sure you have the following:

Part name	Part number
G5478A Microplate Exchanger (single- or dual-stage)	G5508-60011
G5473A Power supply	G5508-60008
Base mounting screws:	
M5 socket-head cap screws (4)	G5550-02377
M6 socket-head cap screws (2)	G5550-02412
Power supply mounting:	
Pan-head Phillips machine screws (4)	G5550-09078
M5 split lock washer (4)	G5550-02453
M5 flat washer (4)	G5550-02439
Power cable	Varies by country
Microplate Exchanger base power cable	G5508-60005
Serial connection:	
Ethernet cable	G5550-22527
Serial adaptor	G5550-21721
<i>Microplate Exchanger User Guide</i>	G5478-90004

If you already have a Microplate Exchanger and you ordered the kit to upgrade the device to a Dual-Stage Microplate Exchanger, make sure you have the following:

Part name	Part number
G5478A Microplate Exchanger (single-stage)	G5508-60011
Microplate Exchanger base power cable	G5508-60005
Serial connection:	
Ethernet cable	G5550-22527
Serial adaptor	G5550-21721
Support posts (4)	5023-1665
M5 socket-head cap screws (8) for the posts	G5550-08530

About installation and setup

Before installing the Microplate Exchanger, make sure the installation site meets the requirements specified in the *Microplate Exchanger User Guide*.

For installation and setup instructions, see the *Microplate Exchanger User Guide*.

About storage and transport

If you need to remove the Microplate Exchanger from the attachment surface and package it for storage or transport, see the *Microplate Exchanger User Guide* for instructions.

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 control software.
- *PDF files*. Product manuals are provided as PDF files on the automation software CD.
- *Agilent Technologies website*. You can search the Automation Solutions product knowledge base or download the latest version of any PDF file 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
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