

# G5581A Microplate Labeler

## Unpacking Guide

This guide describes how to unpack the Microplate Labeler. For installation and setup instructions, see the [G5581A Microplate Labeler User Guide](#).

### Unpacking the shipping box

#### WARNING

#### Before you begin

The Microplate Labeler weighs 18.7 kg (41.3 lb). Do not attempt to unpack or move the Microplate Labeler without assistance. Always use proper lifting techniques to avoid injury.

#### CAUTION

Make sure the shipping box is oriented with the top side up to prevent potential equipment damage.

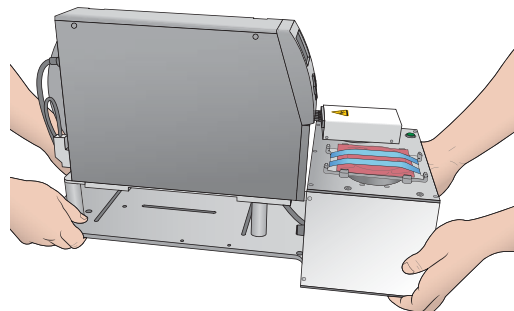
- **Personnel requirements.** Make sure two people are available to lift the Microplate Labeler from the shipping box.
- **Site specifications.** Ensure the installation site meets the required site specifications. For details, see the [G5581A Microplate Labeler User Guide](#).

#### IMPORTANT

Save the packaging materials in case you must move or ship the Microplate Labeler.

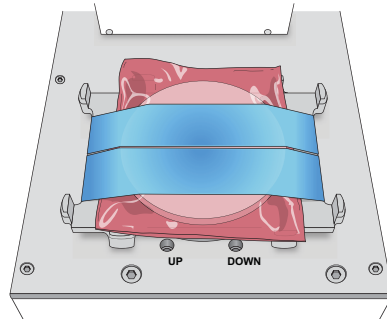
#### To unpack the Microplate Labeler:

- 1 Open the top flaps of the shipping box, and retrieve the Important Documents envelope. Set it aside temporarily.
- 2 Remove the foam packing from the shipping box.
- 3 Carefully unwrap the plastic sheet that covers the Microplate Labeler in the shipping box.
- 4 With a person positioned on each side, reach down inside the shipping box until you can grasp the baseplate of the Microplate Labeler. While firmly holding the baseplate, as the following figure shows, lift the Microplate Labeler onto a lab bench.



## Inspecting the contents

- 5 Remove the round spacer that is taped to the plate stage. Set the spacer aside for potential future use.



### IMPORTANT

To determine if the plate spacer should be installed, you must first complete the device installation and setup described in the *G5581A Microplate Labeler User Guide*. If you determine that the plate spacer is required, contact Agilent Technical Support for installation.

### Inspecting the contents

Inspect all items for completeness and potential shipping damage. If an item is defective on arrival, contact Agilent as soon as possible. For contact information, go to <https://www.agilent.com/en/contact-us/page>.

Verify that the following items are included with your shipment. The items may be packaged in a separate box:

- Air connection kit (non-metric or metric), including tubing and fittings
- Microplate Labeler Media Kit, which includes a roll of labels, thermal transfer ribbon, and a pen for cleaning the printhead.
- Serial and Ethernet communications cables
- Hex wrench (2.5 mm)
- Power cord appropriate for your country

### Accessing user documentation

You can find the automation product user documents in the VWorks Knowledge Base, which is available in the following ways:

- Help menu in the VWorks software
- Online at [www.agilent.com](http://www.agilent.com)/[www.agilent.com/chem/askb](http://www.agilent.com/chem/askb)

### Contacting Agilent

Web: <https://www.agilent.com>

Contact page: <https://www.agilent.com/en/contact-us/page>

