

Vertical Pipetting Station Unpacking Guide

Shipping containers

The Vertical Pipetting Station ships in the following containers:

- **Vertical Pipetting Station crate.** Contains the Vertical Pipetting Station device.
- **Peripherals box.** Contains the following packages:
 - Utility kit
 - Pipette head box
 - Computer box

Depending on the configuration ordered, additional packages or items can be included, such as accessories.

Before you start

Verify the following:

- 1 **Utility kit contents.** Verify the kit includes:
 - Power cable and serial communications cable
 - Tubing and quick-disconnect fittings for air connection (1/4-in) and vacuum (3/8-in), if applicable
 - Software CD-ROM
 - *Vertical Pipetting Station User Guide*
- 2 **Site specifications.** Ensure the installation site meets the requirements specified in the *Vertical Pipetting Station User Guide*.
- 3 **Required tools and equipment.** Obtain the following:
 - Large flat-head screwdriver
 - Cart for moving the device
- 4 **Personnel.** Make sure two people are available for lifting the device from the crate.

WARNING The Vertical Pipetting Station weighs 37.195 kg (82 lb). Do not attempt to move the device without assistance. Use proper lifting techniques when lifting the device.

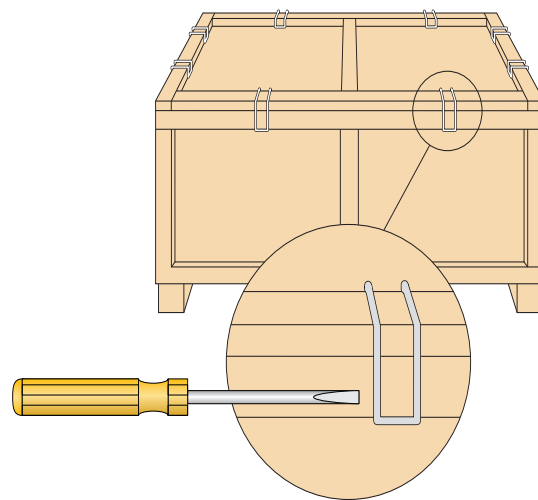
Unpacking the crate

IMPORTANT Save the packaging materials in case you are required to move or ship the Vertical Pipetting Station.

To unpack the Vertical Pipetting Station crate:

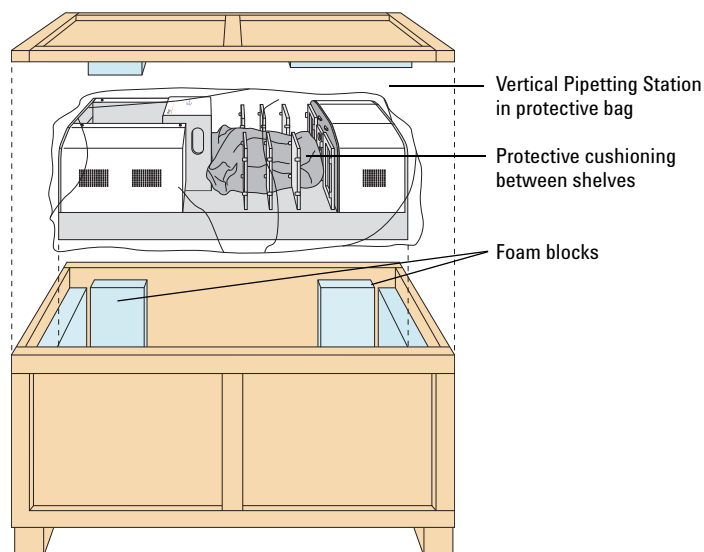
- 1 Use a large flat-head screwdriver to pry off the crate-locking clips, and remove the top of the crate.

WARNING Use care when removing the crate clips. The clips are spring loaded and can pop off when loosened.



CAUTION Do not lift the Vertical Pipetting Station from the shelves, head mount, or any other moving parts. Doing so can damage the device.

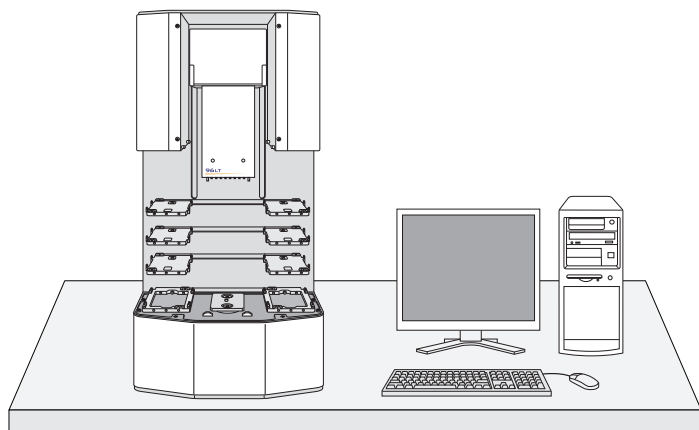
- 2 With a person on either side of the crate, reach between the foam blocks until you can grasp underneath the Vertical Pipetting Station.



- 3 While firmly supporting underneath the Vertical Pipetting Station, lift it up and out of the crate.
- 4 Place the Vertical Pipetting Station in the upright position on the lab bench or table.
- 5 Remove the protective plastic bag covering the Vertical Pipetting Station, and remove the protective cushioning from between the shelves.

About installation and setup

For installation and setup requirements and instructions, see the *Vertical Pipetting Station User Guide*.



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

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 12081443513
euroservice.automation@agilent.com

Documentation feedback:
documentation.automation@agilent.com

Web: www.agilent.com