

Double-Up Gel Drying Rack

Safety

Caution/Warning



The Double-up Gel Dryer Rack is meant to accommodate two gel dryers at once. If only one dryer is placed on the Double-up Rack, it must be placed on the upper shelf. If the gel dryer is placed on the bottom shelf, the Double-up Rack will become unbalanced and tip over, creating a hazardous condition.

Unpacking

The Double-up Gel Dryer Rack is shipped fully assembled. To protect the handle for the pull-out bottom shelf, it is shipped attached to the inside. Remove the two screws attaching the handle using a Philips head screw-driver. Reattach the handle to the front of the shelf.

Set-up

The Double-up Gel Dryer Rack is fully plumbed for vacuum. To set up the Double-up Rack:

1. Place one gel dryer on each shelf.
2. Keeping the tubing in your system to a minimum, attach the vacuum line labeled “UPPER” to the gel dryer on the upper shelf.
3. Attach the vacuum line labeled “LOWER” to the gel dryer on the lower shelf. Be sure to leave enough length of tubing to allow the shelf to be fully extended.
4. Attach the vacuum line labeled “PUMP” to your vacuum source.

Instructions

To operate the gel dryer, follow the instructions in the dryer’s user guide. Check that the valve controls on the front of the rack are in the closed position. Turn on the vacuum source, then open the valves to the gel dryers allowing one gel dryer to seal and then the other.