

3 L and 12 L Harvestainer BPC System

Quick Start Guide

This guide provides a brief overview of 3 and 12 L Thermo Scientific™ Harvestainer™ BPC systems. It is intended to be used by individuals who are already familiar with the system. Prior to using this product, it is important to read and understand the user's guide for detailed warnings, safety, instructions, specifications, and general information.

Warnings and safety

For full details and descriptions of warnings, see the user's guide.

- **Do not** exceed 0.5 psi static BPC pressure
- **Do not** exceed 5 psi inlet pressure
- Static electricity may build up
- Use caution when handling
- Keep sharps away from BPCs

Unpacking information

1. Inspect the packaging for any damage. Call your sales representative if you discover any damage or issues.
2. Open the box carefully to avoid cutting the product.
3. Remove the double polybags, keeping the tray and BPC together.

Preparation and resources

Separate cells and microcarrier beads, as defined by your processes. Then determine the next step in the process. Obtain the following:

- 2 Peristaltic pumps
- Sterile tubing welder and sealer for C-Flex™ 374 tubing 0.95 x 1.59 cm (3/8 x 5/8 in.)—see the specifications.

Process steps

1. Place the Harvestainer BPC on a flat surface, free of sharps (Figure 1).

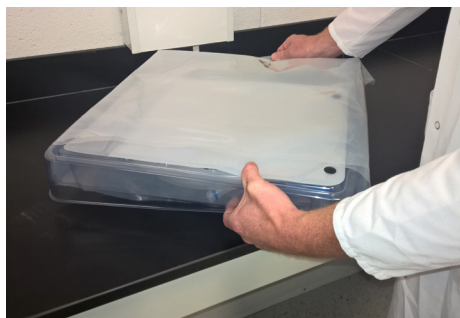


Figure 1. Placing the Harvestainer.

2. Lift the top tray. Place rods into the slots and press firmly into the top tray (Figure 2).

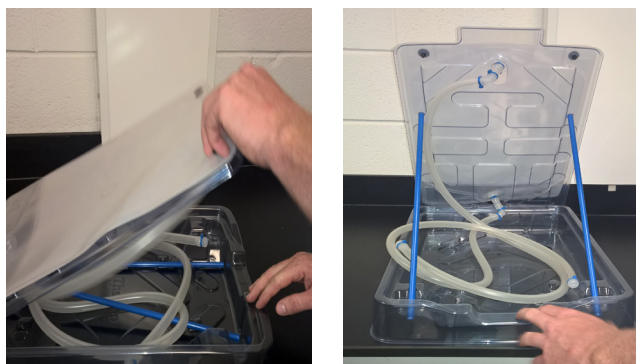


Figure 2. Lifting the tray and placing rods into the slots.

- Remove the polybag from around the tubing lines and extend the tubing to prevent kinking (Figure 3).



Figure 3. Extending the tubing.

- Using the sterile tubing welder, connect the drain line (bottom) to the next step of your process and the transfer line (top) to the bioreactor. Then place the tubing into the pumps (Figure 4).



Figure 4. Connecting the lines and placing tubing in a pump.

- Ensure lines are open. Begin pumping on both lines at 2.5 L/min for 3 L and 3.3 L/min for 12 L.
- When finished, seal off both lines with clamps and/or a tubing sealer.

Complete system setup

Figures 5 and 6 show the complete transfer setup and drain setup, respectively.



Figure 5. Transfer setup, with bioreactor connected to the Harvestainer BPC with pump.



Figure 6. Drain setup, with the Harvestainer BPC connected to storage BPC with pump and scale.

Disposal

Dispose of the entire system, including the single-use support tray, in an appropriate waste container or incinerator.

Specifications

See the table below for an overview of the Harvestainer specifications. See the user's guide for full details and descriptions.

Parameter	3 L	12 L
C-Flex 374 tubing dimensions (inner diameter x outer diameter x length)	0.95 x 1.59 x 122 cm (3/8 x 5/8 x 48 in.)	
Liquid flow for transfer and draining	2.5 L/min	3.3 L/min
Maximum internal BPC pressure rating	0.5 psi	
Microcarrier bead capture capacity	3 L	12 L

Find out more at [thermofisher.com](https://www.thermofisher.com)