

Quick Guide

Dispensing Media and Filling Sample Bags on the ReadyStream® System

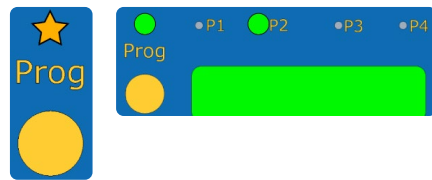
Choose one of 3 possible dispensing parameters for media:

- Use a preset program
- Scan a test data matrix code
- Manually customize the parameters

Preset Program

Press the [PROG] button to move from one preset to another. The LED corresponding to the selected preset will turn green and the respective parameters are displayed.

The system is ready to dispense media.



Manually customize the parameters

To modify the volume, temperature or concentration, the user can select the parameter to be modified by pressing either [ml], [°C] or [X:1].

The system is ready to dispense media. Any customized program can now be stored as a new preset



Dispensing media

Position a container or the test bag in the holder on the dispensing unit. The test bag can be maintained by 2 sticky tack dots (not provided) positioned at the top of each internal face of the holder.

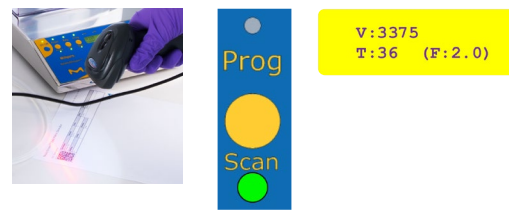
Press [MEDIA] to dispense media according to the displayed parameters of your sample.



Scan a test data matrix code

Scan a test data matrix, the corresponding "Scan" LED turns green and the respective parameters are displayed.

The system is ready to dispense media.



Filling Sample Bags

Select either the standard ReadyStream® bag holder for single samples or the large ReadyStream® bag holder for pooling size samples, and add to the ReadyStream® dispensing unit. The height of the dispensing holder can be adjusted on the stand to accommodate for the bag holders and bag sizes. Standard positions for the standard and large bag holders are marked by yellow tags.



Installation of ReadyStream® consumables

Installation of the ReadyStream® filter set

- 1** Remove the blue cap and connect the filter to the outlet of the media preparation unit



- 2** Install the tubing on the two pins



- 3** Using the barcode reader, scan the data matrix code on the label



Installation of the ReadyStream® media bag

- 1** Place a bag on the media unit. Extend the tray for the 10 L bag



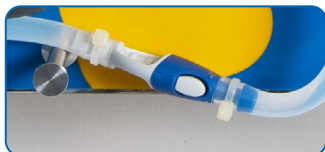
- 2** Install the gray dispensing connector on the media dispensing unit



- 3** Remove the orange cap from the filter and store it in a sterile container



- 4** Connect the blue connector of the media bag to the grey connector of the filter set



- 5** Insert the stopcock valve in the actuator



- 6** Position the tubing in the guiding pins



- 7** Position the tubing through the peristaltic pump. Do not close the pump



- 8** Using the barcode reader, scan the data matrix code on the media bag



To place an order or receive technical assistance

Order/Customer Service:
[SigmaAldrich.com/order](https://www.sigmaaldrich.com/order)

Technical Service:
[SigmaAldrich.com/techservice](https://www.sigmaaldrich.com/techservice)

MilliporeSigma
400 Summit Drive
Burlington, MA 01803

