

# QUICK START GUIDE

## STEMprep™ Tissue Dissociator

For detailed operating instructions, please refer to the Technical Manual for STEMprep™ Tissue Dissociator (Document #10000030598) available at [www.STEMCELL.com](http://www.STEMCELL.com).



### 1 Start Up STEMprep™ Tissue Dissociator

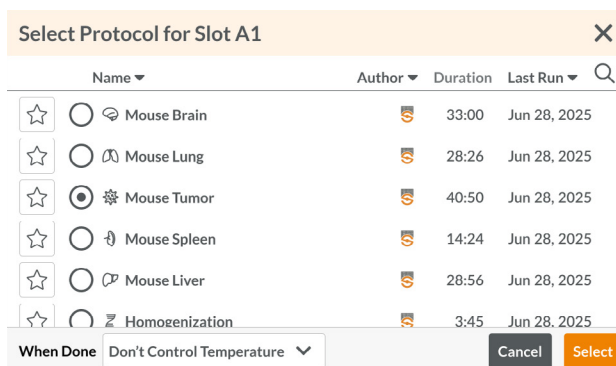
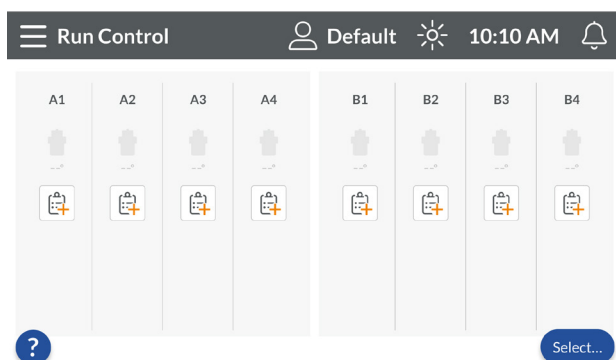
Start up STEMprep™ Tissue Dissociator using the **power switch** located on the back of the instrument. The touch screen on the Primary Unit will start up and the unit label will appear on each connected unit.



### 2 Select Desired Protocol

After startup, the **Run Control** screen will display a representation of the slots for each installed unit. In the example below, 8 slots are available for two installed STEMprep™ units: unit A (slots A1 - A4) and unit B (slots B1 - B4).

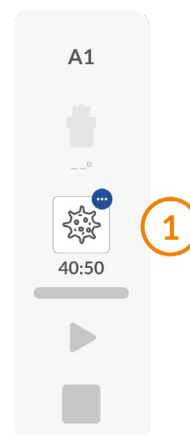
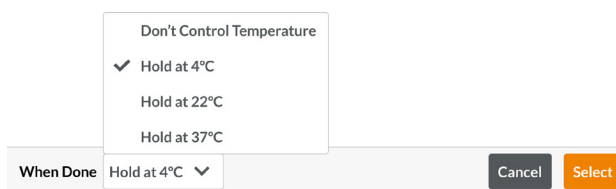
- Choose the slot you would like to use by tapping the **Select Protocol** button .
- A list of available protocols will be displayed. Tap the desired protocol name to select it. The selected protocol is indicated with a filled-in radio button .



iii) **Optional:** Tap the **When Done** dropdown menu to set a post-protocol hold temperature.

iv) Tap **Select** to apply the selected protocol and hold temperature to the slot.

v) The tissue type of the selected protocol will now be displayed for that slot ①. A blue circle indicates the application of a post-protocol temperature hold.



### 3 Prepare the STEMprep™ Sample Tube

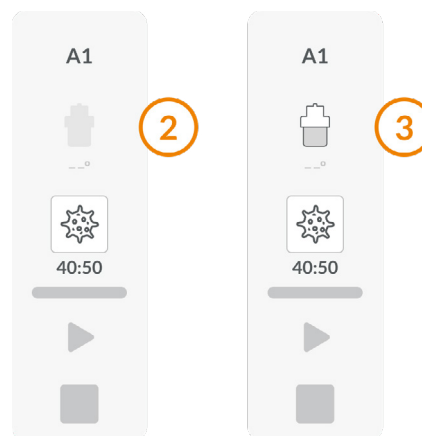
i) Add the tissue sample to a STEMprep™ Sample Tube with the applicable STEMprep™ Tissue Dissociation Kit Cocktail.



### 4 Load the STEMprep™ Sample Tube



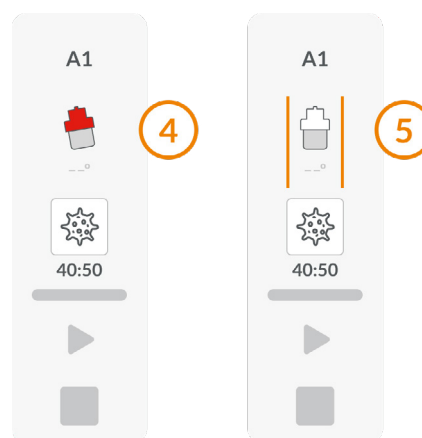
- i) Initially, the **Run Control** screen will display a dimmed **sample tube** icon ② for the empty slot (to indicate that no sample tube is present). Insert a STEMprep™ Sample Tube in the slot; the icon will change to indicate a sample tube is present ③.



- ii) If the tube is not properly inserted, the **sample tube** icon will indicate that it is **misaligned** ④.

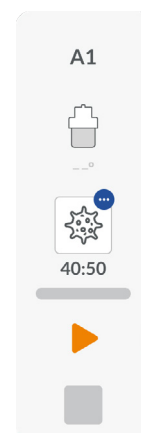
**Note:** When the shield is lowered to its lowest position, the instrument checks whether the spindle is making proper contact with the STEMprep™ Sample Tube. If proper engagement is not detected, the spindle will automatically move back and forth in an attempt to align correctly. If engagement still cannot be achieved, the spindle will stop moving. In this case, manually raise the shield and gently rotate the spindle by a few degrees before trying again.



The **shield raised** indicator ⑤ is displayed while the shield for that slot is lifted, whether or not a sample tube is present.

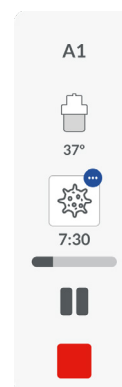


## 5 Start the Protocol

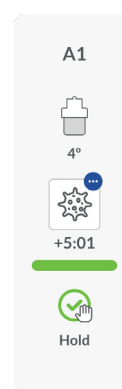
- i) When the sample tube is inserted correctly and the protocol is selected, the **Start** button ► is enabled. Tap the Start button to begin running the protocol.



- ii) Once the protocol is running, the timer will count down and the progress bar will visually display the completion progress. To pause or stop the run, tap the **Pause** button  or the **Stop** button  , respectively.



- iii) Once the protocol run is complete, the timer will count up from the time of completion.



## 6 Unload the STEMprep™ Sample Tube

Remove the sample tube. The sample is now ready for downstream processing.

For detailed instructions on processing various tissues, including brain, lung, tumor, and more, refer to the Product Information Sheet (PIS) for the applicable STEMprep™ Tissue Dissociation Kit. PISs are available at [www.STEMCELL.com](http://www.STEMCELL.com); you may also contact us to request a copy.



## 7 Shut Down STEMprep™

Shut down the instrument using the power switch located at the back of the unit.

Copyright © 2025 by STEMCELL Technologies Inc. All rights reserved including graphics and images. STEMCELL Technologies & Design, STEMCELL Shield Design, Scientists Helping Scientists, and STEMprep are trademarks of STEMCELL Technologies Canada Inc. All other trademarks are the property of their respective holders. While STEMCELL has made all reasonable efforts to ensure that the information provided by STEMCELL and its suppliers is correct, it makes no warranties or representations as to the accuracy or completeness of such information.

UNLESS OTHERWISE STATED, PRODUCTS ARE FOR RESEARCH USE ONLY AND NOT INTENDED FOR HUMAN OR ANIMAL DIAGNOSTIC OR THERAPEUTIC USES. FOR PRODUCT-SPECIFIC COMPLIANCE AND INTENDED USE INFORMATION, REFER TO THE PRODUCT INFORMATION SHEET. GENERAL INFORMATION ON QUALITY AT STEMCELL MAY BE FOUND AT [WWW.STEMCELL.COM/COMPLIANCE](http://WWW.STEMCELL.COM/COMPLIANCE).