

PowerEase™ Touch 120W Power Supply

Cat. no. PS0120 (115 VAC), PS0121 (230 VAC), PS0122 (230 VAC Australia), PS0123 (230 VAC China)

Publication No. MAN0019261

Rev. B.0

Use the following instructions to operate the PowerEase™ Touch 120W Power Supply in constant voltage mode. For more information and detailed instructions, see the power supply manual available at thermofisher.com.

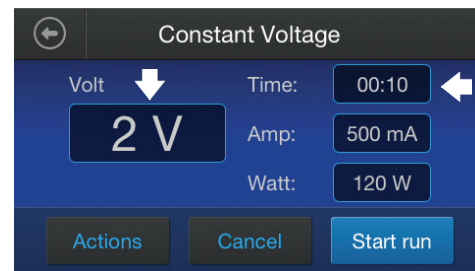
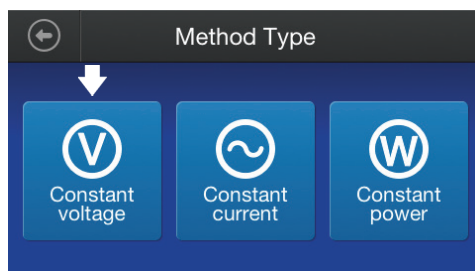
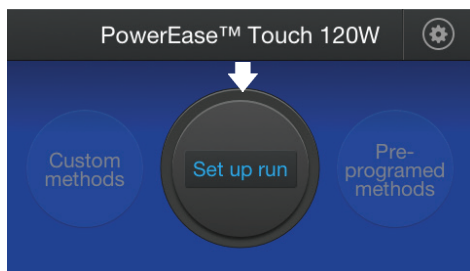
Set up the power supply

1. Check the label located by the AC inlet of the power supply to ensure that the unit is compatible with locally provided voltage.
2. Place the power supply on a level laboratory bench in an area with proper ventilation. Position the unit such that the **On/Off** switch and the AC inlet at the rear of the power supply are easily accessible.
3. Ensure the **On/Off** switch located at the rear of the power supply is in the **Off** position.
4. Attach the power cord to the AC inlet of the power supply.
5. Plug the power supply into an appropriate electrical outlet (use properly grounded AC outlets and power cords).
6. Connect the leads from an electrophoresis unit to the power supply by inserting the positive lead (+) into the red output jack and the negative lead (-) into the black output jack.
7. Turn the AC power switch at the back of the power supply **On**.

Note: When the power supply is initially switched on, **Set up run**, **Custom methods**, and **Pre-programmed methods** are displayed as options on the home screen. After a saved or pre-programmed method is used, **Last run** is displayed as an option in the center dial.

Operate the power supply (constant voltage)

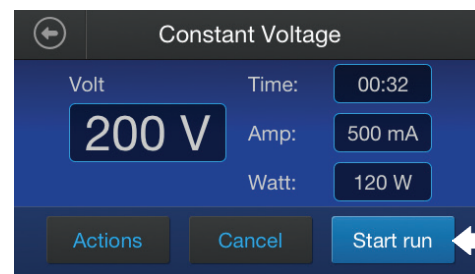
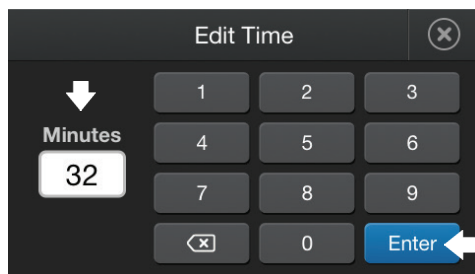
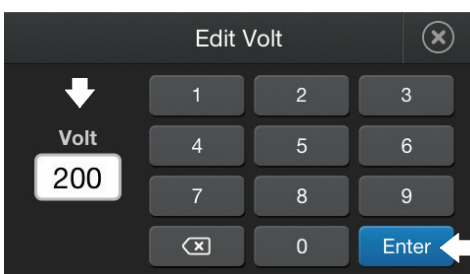
The PowerEase™ Touch 120W Power Supply is a medium throughput power supply that uses a touch-screen user interface. Constant Voltage Mode is recommended when using the instrument with protein gels or blots. For instructions using other operating modes see the power supply manual available at thermofisher.com.



1. Turn on the power supply, then select **Set up run**.

2. Select the **Constant voltage** operating mode for **Method Type**.

3. Select the data fields for the run parameters to be changed from the default values. ^[1]



4. Set the operating mode parameter (**Constant voltage**) by entering the value in the **Volt** field, then select **Enter**.

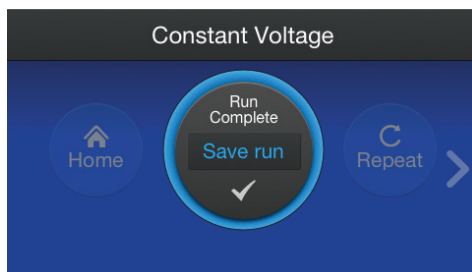
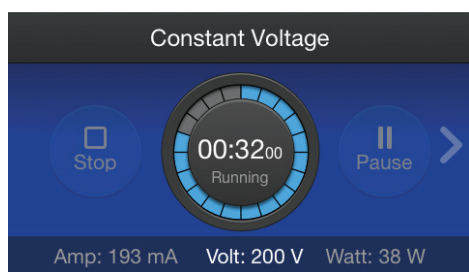
5. Set the time parameter by entering the value in the **Time** field, then select **Enter**.

6. Select **Start run** to initiate the method.

[1] When editing constant mode parameters, it is recommended that only the operating mode and the time parameter are changed, leaving other parameters at the default power supply maximums in order to prevent unwanted limitations in power delivery during the run.

Continued on reverse side.

Operate the power supply (constant voltage), continued



7. The run screen displays the run parameters in real time during the run. The center dial displays a countdown timer for the method.

Select the arrow (➤) to display run settings.

8. Completion of the run when the timer reaches 00:00:00 is indicated by an audible beeping alarm and a display of **Run Complete** ✓ in the center dial.

Select **Home** to return to the home screen.

Select **Repeat** to re-run the same method.

Select **Save run** to save the method settings for future use (this option is not available if the method has already been saved).

Limited Product Warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale at www.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have any questions, please contact Life Technologies at www.thermofisher.com/support.



Life Technologies | Carlsbad, CA 92008 USA | Toll Free in USA 1.800.955.6288

For descriptions of symbols on product labels or product documents, go to thermofisher.com/symbols-definition.

The information in this guide is subject to change without notice.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Important licensing information: This product may be covered by one or more Limited Use Label Licenses. By use of this product, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2021 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

For support visit thermofisher.com/support | thermofisher.com/askaquestion

thermofisher.com

04 June 2021

ThermoFisher
SCIENTIFIC