

QUICK REFERENCE

PowerEase® 90W Power Supply

Catalog Numbers PS0090, PS0091, PS0092

Publication Number MAN0009567

Rev. 1.0

Use the following instructions to install and operate the PowerEase® 90W Power Supply. For more information and detailed instructions, refer to the *PowerEase® 90W Power Supply User Guide*, available at www.lifetechnologies.com/manuals.

Note: For safety guidelines, refer to the "Safety" appendix in the *PowerEase® 90W Power Supply User Guide*.

Install the PowerEase® 90W Power Supply

1. Check the label located near the AC inlet to help ensure that the unit is compatible with locally provided voltage.
2. Place the PowerEase® 90W Power Supply on a level laboratory bench, in an area with proper ventilation.
3. *For your safety:* Position the unit properly such that the **On/Off** switch and the AC inlet located on the rear of the unit are easily accessible.
4. Ensure the AC power switch is in the **Off** position and attach the power cord to the AC inlet of the power supply.
Note: Use only properly grounded AC outlets and power cords.
5. Plug the power supply into the appropriate electrical outlet.
6. Connect the leads from an electrophoresis unit to the power supply by inserting the positive lead (+) into the positive (red) output jack, and the negative lead (-) into the negative (black) output jack.

PowerEase® 90W Power Supply key pad

Front view of the PowerEase® 90W Power Supply.



Press...	To...
^ (Up Arrow)	Move the cursor up and increase values when settings are flashing.
v (Down Arrow)	Move the cursor down and decrease values when settings are flashing.
Mode	Cycle between voltage, current, and time settings, allowing values to be changed.
Constant	Toggle between constant voltage or current operation.
Start/Pause	Start, pause, and resume power supply operation without resetting the timer.
Stop	Stop power supply operation <i>and</i> reset the timer.

Quick operation instructions

Use the following instructions to operate the PowerEase® 90W Power Supply.

1. Turn the AC power switch at the back of the power supply to the **On** position.
 - The screen displays the last set **V** (voltage) in large bold type on the left side of the display and **TIME** and **AMPs** in small type on the right side of the display.
2. Press **Constant** to toggle between Constant Voltage Mode and Constant Current Mode.
 - The screen displays the selected operating mode in large bold type on the left of the display.
3. Press **Mode** to edit the parameters of the operating mode: Volts, Hours, Minutes, and Amps.
4. While a value is flashing, you may use the directional arrow buttons (v/^) to change the values.
 - The values stop flashing and the instrument exits editing mode 5 seconds after the last key has been pressed.
 - If you set **Time** to **00:00**, the power supply will run continuously.
5. Press **Start/Pause** to start electrophoresis.
 - The red light on the display will illuminate while power is applied to the load.
 - Electrophoresis stops when the timer reaches **00:00** if a time has been set.
 - When electrophoresis stops, the power supply *beeps* and the screen flashes **TIME UP**.
6. Press **Stop** to silence the *beep*.

For Research Use Only. Not for use in diagnostic procedures.



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 found on Life Technologies' website at www.lifetechnologies.com/termsandconditions. If you have any questions, please contact Life Technologies at www.lifetechnologies.com/support.

For Research Use Only. Not for use in diagnostic procedures.

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

Limited Use Label License: Research Use Only: The purchase of this product conveys to the purchaser the limited, non-transferable right to use the purchased amount of the product only to perform internal research for the sole benefit of the purchaser. No right to resell this product or any of its components is conveyed expressly, by implication, or by estoppel. This product is for internal research purposes only and is not for use in commercial applications of any kind, including, without limitation, quality control and commercial services such as reporting the results of purchaser's activities for a fee or other form of consideration. For information on obtaining additional rights, please contact outlicensing@lifetech.com or Out Licensing, Life Technologies, 5791 Van Allen Way, Carlsbad, California 92008.

DISCLAIMER: LIFE TECHNOLOGIES CORPORATION AND/OR ITS AFFILIATE(S) DISCLAIM ALL WARRANTIES WITH RESPECT TO THIS DOCUMENT, EXPRESSED OR IMPLIED, INCLUDING BUT NOT LIMITED TO THOSE OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, OR NON-INFRINGEMENT. TO THE EXTENT ALLOWED BY LAW, IN NO EVENT SHALL LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) BE LIABLE, WHETHER IN CONTRACT, TORT, WARRANTY, OR UNDER ANY STATUTE OR ON ANY OTHER BASIS FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING BUT NOT LIMITED TO THE USE THEREOF.

©2013 Life Technologies Corporation. All rights reserved. The trademarks mentioned herein are the property of Life Technologies Corporation and/or its affiliates or their respective owners.

Headquarters

5791 Van Allen Way | Carlsbad, CA 92008 USA | Phone +1 760 603 7200 | Toll Free in USA 800 955 6288

For support visit lifetechnologies.com/support

lifetechnologies.com

17 September 2013

