

PowerEase® 300W Power Supply

Catalog Numbers PS0300, PS0301, PS0302

Publication Number MAN0008921

Rev. 1.0

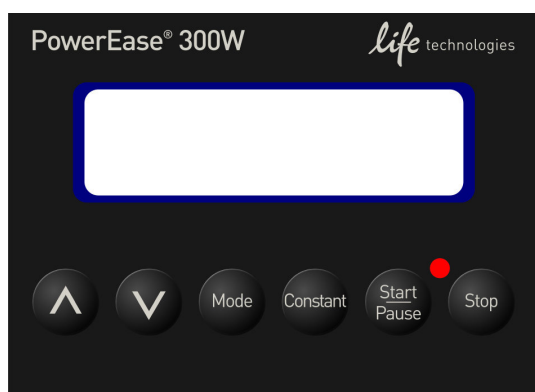
Use the following instructions to operate the PowerEase® 300W Power Supply. For more information and detailed instructions, refer to the PowerEase® 300W Power Supply manual, available at www.lifetechnologies.com/manuals.

Before starting

1. Ensure the AC power switch located on the back of the power supply is in the **Off** position.
2. Attach the power cord to the AC inlet of the power supply.
Note: Use only properly grounded AC outlets and power cords.
3. Plug the power supply into the appropriate electrical outlet.
4. 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.

PowerEase® 300W Power Supply key pad

Front view of the PowerEase® 300W Power Supply.



Press...	To...
^ (Up Arrow)	Move the cursor up and increase values when settings are flashing.
∨ (Down Arrow)	Move the cursor down and decrease values when settings are flashing.
MODE	Select Constant Mode or Program Mode . Note: The MODE key cycles between voltage, current, power, and time settings, allowing values to be changed.
CONSTANT	Toggle between constant voltage, current, or power 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.

Constant mode quick operation instructions

Use the following instructions to operate the PowerEase® 300W Power Supply in Constant Mode. For Program Mode instructions, refer to the PowerEase® 300W Power Supply manual.

Note: The PowerEase® 300W is a high throughput power supply. For optimal performance, Life Technologies recommends using Constant Voltage Mode (see step 3, below) when using the instrument with 1 or 2 protein gels or blots.

1. Turn the AC power switch at the back of the power supply to the **On** position.
 - The PowerEase® 300W Power Supply screen displays **Constant Mode** and **Program Mode**.
 - **Constant Mode** will be displayed in yellow letters surrounded by a black box.

Note: You may use the directional arrow buttons (∨/∧) to toggle between **Constant Mode** and **Program Mode**, if needed.
2. Press **MODE** to select constant mode operation.
 - The Power Ease 300W Power Supply defaults to Constant Voltage Mode operation.
 - The screen displays the last set **V** (voltage) in large bold type on the left side of the display and **TIME**, **AMPS**, and **WAT** (watts) in small type on the right side of the display.
3. Press **CONSTANT** to toggle between the operating modes: Constant Voltage Mode, Constant Current Mode, and Constant Power Mode.
 - The screen displays the selected operating mode in large bold type on the left of the display.
4. Press **MODE** to edit the parameters of the operating mode: Volts, Hours, Minutes, Amps and Watts.
5. While a value is flashing, you may use the directional arrow buttons (∨/∧) 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.
6. 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**.
7. Press **STOP** to silence the *beep* and return to the Constant Operating Mode display.

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.

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.

©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

16 September 2013

