

## ELECTROPHORESIS

# PowerPac Universal Power Supply Method Mode Operation

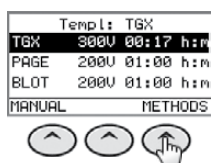
### Quick Guide

This quick guide describes how to program and perform a run in Method mode. Method mode operation allows users to create and run single-step and multistep methods. In method mode, up to nine methods, each with up to nine steps, can be programmed, stored, or edited. Refer to the PowerPac Universal Power Supply Instruction Manual (#4006223) for complete instructions.

## Create and Run New Method

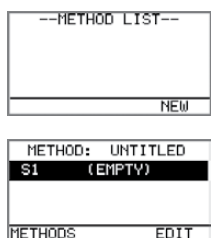
### 1 Power on instrument.

### 2 Select the programming mode.



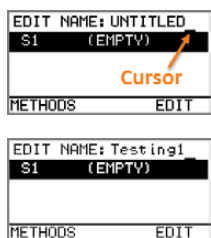
On the Home screen, press **METHODS** to display the Method List screen.

### 3 Create a new method.



To create a new method, press **NEW**.

### 4 Name the new method.

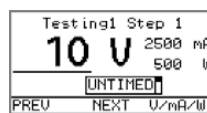


Press the up arrow key and a blinking cursor appears.

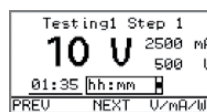
Delete existing name (UNTITLED) by pressing **CE** (on numeric soft keypad) and enter the new method name.

Press the down arrow key to accept the method name.

### 5 Edit the first step in the method.

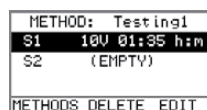


Press **EDIT** (on numeric soft keypad) and select volts (V), milliamps (mA), or watts (W).



Adjust the run parameters:

- V: 10–2,500
- mA: 10–2,500
- W: 1–500
- V hours (timer): 1–99,999
- hh:mm (timer): 00:01–99:59

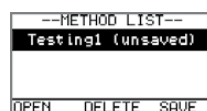


Press **EDIT** to navigate back to the Method screen to add steps if necessary.

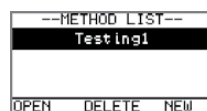
### 6 Save the method.



On the Method screen, press **METHODS** to access the Method List screen.



Press **SAVE** to save the method.



### 7 Start the run.

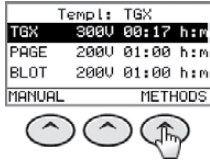


Press the run key to start the run.

# Edit (or View) and Run a Saved Method

**1 Power on instrument.**

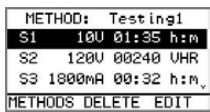
**2 Select the programming mode.**



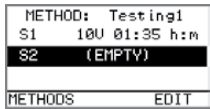
On the Home screen, press **METHODS** to display the Method List screen.



Select a method to edit.



Press the up or down arrow key to select the step you want to edit, then press **EDIT** to view the step parameters.



Add steps to the method by following the instructions in step 5 of the Create and Run New Method section.

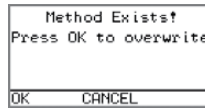
**3 Save the method.**



On the Method screen, press **METHODS** to access the Method List screen.



Press **SAVE** to save the method.



Press **OK** to save changes.



**4 Start the run.**



Press the run key to start the run.

Visit [bio-rad.com/PowerSupply](http://bio-rad.com/PowerSupply) for more information.

Bio-Rad is a trademark of Bio-Rad Laboratories, Inc. in certain jurisdictions. All trademarks used herein are the property of their respective owner.



**Bio-Rad  
 Laboratories, Inc.**

Life Science  
 Group

**Web site** [bio-rad.com](http://bio-rad.com) **USA** 1 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 43 01 877 89019 **Belgium** 32 03 710 53 00 **Brazil** 55 11 3065 7550  
**Canada** 1 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 36 01 459 6192 **Denmark** 45 04 452 10 00 **Finland** 35 08 980 422 00 **France** 33 01 479 593 00  
**Germany** 49 089 3188 4393 **Hong Kong** 852 2789 3300 **Hungary** 36 01 459 6190 **India** 91 124 4029300 **Israel** 972 03 963 6050 **Italy** 39 02 49486600  
**Japan** 81 3 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 31 0 318 540 666 **New Zealand** 64 9 415 2280 **Norway** 47 0 233 841 30  
**Poland** 36 01 459 6191 **Portugal** 351 21 4727717 **Russia** 7 495 721 14 04 **Singapore** 65 6415 3188 **South Africa** 36 01 459 6193 **Spain** 34 091 49 06 580  
**Sweden** 46 08 555 127 00 **Switzerland** 41 0617 17 9555 **Taiwan** 886 2 2578 7189 **Thailand** 66 2 651 8311 **United Arab Emirates** 971 4 8187300  
**United Kingdom** 44 01923 47 1301

