

## Opening and Closing the Heated Lid

### Opening the Lid

1. Turn the knob counterclockwise to raise the inner lid.
2. Raise the handle up to the vertical position.



### Closing the Lid

1. Using the handle, push the lid down.
2. Lower the handle to the horizontal position until the lid latches securely (see inset).
3. Turn the knob clockwise until the inner lid contacts the vessels, then add one-half turn for sealing tape or one-quarter turn for caps.



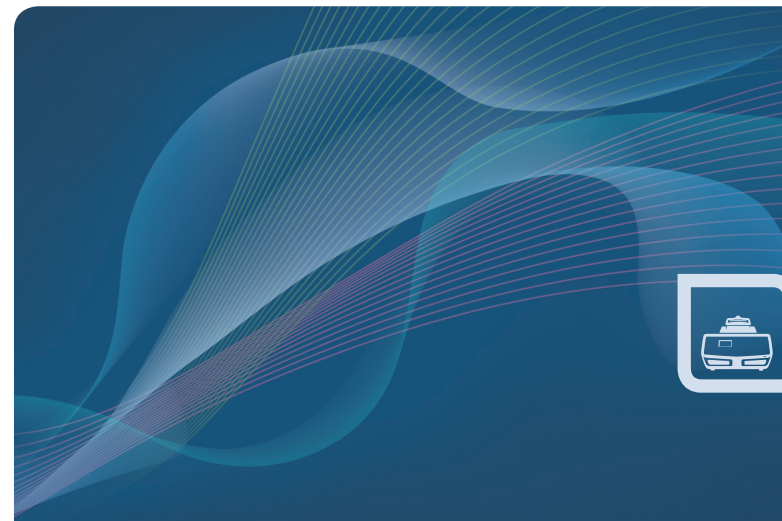
Purchase of this instrument conveys a limited non-transferable immunity from suit for the purchaser's own internal research and development and for use in human in vitro diagnostics and all other applied fields under U.S. Patent Number 5,475,610 (Claims 1, 44, 158, 160-163, and 167 only), or corresponding claims in its non-U.S. counterpart, owned by Applied Biosystems. No right is conveyed expressly, by implication, or by estoppel under any other patent claim, such as claims to apparatus, reagents, kits, or methods such as 5' nuclease methods. Further information on purchasing licenses may be obtained by contacting the Director of Licensing, Applied Biosystems, 850 Lincoln Centre Drive, Foster City, California 94404, USA.

This product is covered by one or more of the following U.S. patents or their foreign counterparts owned by Eppendorf AG: U.S. Patent Numbers 6,767,512 and 7,074,367.


**BIO-RAD**
**Bio-Rad  
Laboratories, Inc.**

 Life Science  
Group

**Web site** [www.bio-rad.com](http://www.bio-rad.com) **USA** 800 424 6723  
**Australia** 61 2 9914 2800 **Austria** 01 877 89 01 **Belgium** 09 385 55 11  
**Brazil** 55 31 3689 6600 **Canada** 905 364 3435 **China** 86 21 6169 8500  
**Czech Republic** 420 241 430 532 **Denmark** 44 52 10 00  
**Finland** 09 804 22 00 **France** 01 47 95 69 65 **Germany** 089 31 884 0  
**Greece** 30 210 777 4396 **Hong Kong** 852 2789 3300  
**Hungary** 36 1 459 6100 **India** 91 124 4029300 **Israel** 03 963 6050  
**Italy** 39 02 216091 **Japan** 03 6361 7000 **Korea** 82 2 3473 4460  
**Malaysia** 60 3 2117 5260 **Mexico** 52 555 488 7670  
**The Netherlands** 0318 540666 **New Zealand** 64 9 415 2280  
**Norway** 23 38 41 30 **Poland** 48 22 331 99 99 **Portugal** 351 21 472 7700  
**Russia** 7 495 721 14 04 **Singapore** 65 6415 3170  
**South Africa** 27 861 246 723 **Spain** 34 91 590 5200  
**Sweden** 08 555 12700 **Switzerland** 061 717 95 55  
**Taiwan** 886 2 2578 7189 **Thailand** 66 2 6518311  
**United Kingdom** 020 8328 2000

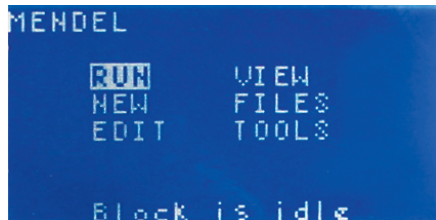
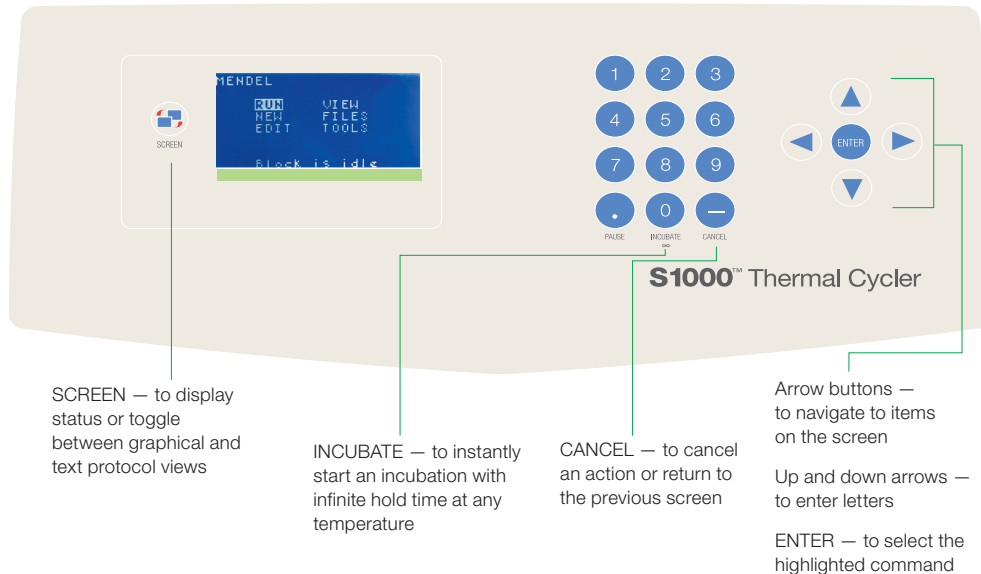


## S1000™ Thermal Cycler Quick Guide

**BIO-RAD**

## Main Menu and Controls

Use the main menu to access the thermal cycler's major functions: RUN a protocol, create a NEW protocol, EDIT a protocol, VIEW a protocol, select FILES, or access TOOLS to change the thermal cycler settings. Use the keys on the instrument's control panel for the functions shown below.



Use the main menu to access major functions

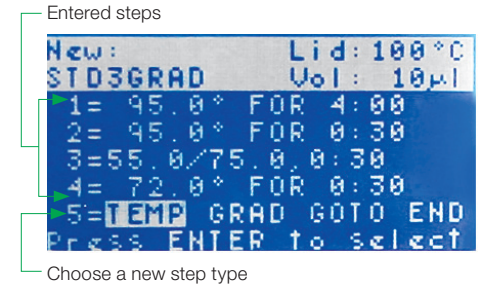
## Running a Protocol

1. Select RUN in the main menu; press ENTER.
2. Select a protocol; press ENTER.
3. Edit the volume, if desired; press ENTER.
4. Select RUN; press ENTER.



## Creating a Protocol

1. Select NEW in the main menu; press ENTER.
2. Enter a name and sample volume; press ENTER.
3. For the initial step, select TEMP for a temperature step or GRAD for a gradient step and enter values; press ENTER.
4. For additional steps, select TEMP, GRAD, or GOTO and enter values; press ENTER.
5. To complete the protocol, select END; press ENTER.
6. Select a folder and save the protocol.



## Editing a Protocol

1. Select EDIT in the main menu; press ENTER.
2. Select a protocol; press ENTER.
3. Edit the sample volume and/or lid temperature, if desired.
4. Use the arrows to highlight and change a value.
5. To insert or delete a step, highlight the step number and select INS or DEL.
6. Select END; press ENTER to finish; changes are saved to the file you selected.

