

www.BenchmarkScientific.com

Benchmark
Scientific 
P.O. Box 709 Edison, NJ 08817 USA
Ph: 908-769-5555 EM: info@BenchmarkScientific.com



Infrared Sterilizer Operations Manual

2011



Benchmark
Scientific 
P.O. Box 709 Edison, NJ 08817 USA
Ph: 908-769-5555 EM: info@BenchmarkScientific.com

Thank you for selecting our BactiZapper infrared sterilizer.

Your BactiZapper has been designed with durability, reliability, and safety in mind. It is your responsibility to install this instrument in conformance with local electrical codes.

This manual contains important operating and safety information. You must carefully read and understand the contents of this manual prior to operating this instrument.

Contents

1.) Introduction.....	1
2.) Technical data.....	2
3.) Operating Instructions.....	3
4.) Precautions and Maintenance	4
5.) Warranty Information.....	5
6.) Included Materials.....	5

5.) Warranty Information

5.1) The BactiZapper Sterilizer unit itself (with the exception of the heater element) is warranted for one year against defects in material and workmanship. The warranted life of the heater element is 3 months.

5.2) Should damage to the instrument occur due to damage, abuse, improper use or improper maintenance, this written warranty is void.

6.) Included Materials

No.	Item	Type	QTY
1	BactiZapper Infrared Sterilizer	B1000	1
2	Operations manual		1
3	Fuse	250V 3A	2

4.) Precautions and Maintenance

4.1) Only insulated loop holders should be used when inserting inoculating loops or needles into the heater element.

4.2) The BactiZapper Sterilizer is not intended for use with scalpels, forceps or any sharp objects. The use of such objects can cause damage and element malfunction and will void the heater element warranty.

4.3) Visual inspection of the heater element should be performed periodically to determine if the heater element core is worn. Inspect for small cracks and residue buildup in both the cool and heated conditions. In the heated condition, cracks can be seen as small, intensely yellow-orange splits. If any defects are noted, the heater element should be replaced before element malfunction occurs.

4.4) The inoculating loops and needles should not be allowed to remain in the heater element for extended periods of time. This will cause the loop holder to become extremely hot and shorten the usable life of the loops and needles.

4.5) Use only the proper fuse for replacement.

4.6) The outer shield of the heater element can reach temperatures as high as 400° F. Do not touch it or introduce any flammable or volatile materials into the general area surrounding the BactiZapper.

4.7) Do not use a two prong adapter to connect the electrical outlet. This instrument must be properly grounded.

1.) Introduction

The BactiZapper Sterilizer is a gasless, flameless sterilizer for use with platinum inoculating loops, needles, glass tube/pipette mouths and various metal & borosilicate glass instruments. Sterilization is accomplished through infrared heat without the dangerous spattering of microorganisms.

Before starting up this unit for the first time, please read the operations manual.

1.1) Structure Description



2.) Technical data

Model	B1000
Power supply	115V / 220 V: 50-60Hz
Power	135W
Highest temperature at core	825°C ± 50°C
Largest possible diameter of item to be sterilized	14mm
Length of heating area	100mm
Heating time	20min
Fuse (US, North America)	250V 3A 5X20mm
Operating temperature	5°C ~ 40°C
Dimensions (W×D×H)	4 x 5 x 7 in. 102mm×130mm×180mm
Net Weight	2.9 lbs. / 1.3kg

3.) Operating Instructions

- 3.1) Remove the BactiZapper from all packaging and place on flat surface.
- 3.2) Plug unit into power outlet. The voltage must correspond to the voltage listed on the product nameplate.

3.3) The red indicator light will come on when the toggle switch is in the “High” position. Twenty minutes after the Bacti-Cinerator Sterilizer is plugged in and the switch is in the “high” position, the optimum sterilizing temperature (825 ° C) will be reached. If the unit is not going to be used for extended periods of time, please set the power switch in “Off” position to conserve the life of the heater element.

Note: To extend the life of the heating element, do not keep the BactiZapper in the high position for more than three hours. If left on for more than three hours, please switch to the “low” position for 30 min. before re-setting the BactiZapper to the high position.

Green Indicator Light – Indicates Low Temp. setting, 500 ° C
Red Indicator Light – Indicates High Temp. setting, 825 ° C

3.4) Gently insert the item to be sterilized into the cylindrical sterilization area. ALWAYS ENSURE THAT THE ITEM IS HELD WITH AN INSULATED MATERIAL. DO NOT DIRECTLY TOUCH THE STERILIZED ITEM! When inserting, avoid scraping the sides of the element to ensure the longevity of your loop and the heater element. The loop should remain inside the heater element area for a minimum of five seconds. Microorganisms are generally destroyed within five seconds. It is not necessary for the needle or loop to be glowing.