



MyTemp™ Digital Mini Incubators

Model H2200-HC and H2200-H Instruction Manual

Benchmark's MyTemp Mini Incubators are micro-processor controlled instruments designed for providing accurate temperatures up to 60°C. With large incubation chambers and removable shelves, the MyTemp incubators are useful for a wide variety of laboratory applications.

I. Product Specifications:

Capacity: 20L

Temperature Range:

H2200-HC: Ambient -15°C to 60°C. (Heating & Cooling Model)

H2200-H 20°C to 60°C (Heat Only Model)

Temperature Accuracy: +/-0.5°C (37°C) **Temperature Uniformity:** +/-1°C (37°C)

Dimensions (Internal): wxdxh 26 x 23.5 x 32.5 cm (10.25 x 9.25 x 12.75")

Dimensions (External): wxdxh 33.5 x 37x 47.5 cm (13.25 x 14.5 x 18.75")

Weight: 6.5 kg (15 lbs)

Warranty: 2 years

Electrical: 100-240VAC, 50/60Hz

II. Product Set-Up:

Place the incubator on a clean, level and stable surface. Plug the power cord into the back of the unit, then plug the electrical cord into a properly grounded outlet. Once plugged in, press the power button on the front display.

III. Product Operation:

To select the desired temperature, press the "Up" or "Down" key. The display begins to flash and now the "Up" and "Down" keys can be used to choose the desired temperature. Once the desired temperature has been selected, the display stops flashing and the chamber begins to heat or cool accordingly.

WARNING: When loading samples, always avoid placing samples directly in front of fan vents. This will minimize airflow and diminish temperature performance.

The incubator is also equipped with an internal light to assist in viewing. To active the internal light, press the "Light" button. To turn off the internal light, press the "Light" button again.



V. Care and Maintenance:

No routine maintenance is required other than to keep the instrument clean. When cleaning, use a damp cloth to wipe down the internal or external parts of the incubator. Avoid the use of solvents as they may attack the product housing. In addition, always ensure that the fan vents are clear of dirt or any other obstructions.

Website: www.BenchmarkScientific.com

Tel: 908-769-5555

Email: info@benchmarkscientific.com

Fax: 908-222-1864