



Mortexer™ Vortex Mixer Model BV1005(E) Instruction Manual

The Mortexer™ is MORE than your standard vortexer mixer. The unique design of the Multi-Head™ (included) features a traditional cup head for general purpose mixing, plus holds up to eight microcentrifuge tubes. Benchmark's Mortexer™ is a quiet, very powerful, vortex mixer designed for vortexing micro tubes of all sizes (0.4ml to 50ml and larger).

I. Product Specifications:

Speed / Control: User settable 200 to 3400 rpm

Modes: ON (continuous), OFF and TOUCH mix

Head: Multi-Head™ for general mixing and up to 8 microtubes

Dimensions (wxdxh): 130x160x170mm, Weight: 3.8kg, 8.4lb

Electrical: 115VAC, 60Hz, 1.5 amps (Model BV1005)

230VAC, 50Hz, 0.75 amps (Model BV1005-E) **CE** compliant

230VAC, 60Hz, 0.75 amps (Model BV1005-E2) **CE** compliant

Operating environment: 4° C to 65° C;

Warranty: 2 years

II. Product Set-Up:

Place the vortex mixer on a clean, flat, stable surface and press down slightly to allow the suction cup feet to adhere. Make sure that the mode switch is in the OFF position and plug the vortex mixer into a properly grounded outlet. Using the SPEED control knob, adjust the mixing speed to the desired setting. Start at the middle of the speed range if you are uncertain of what setting you need. Check that the cup head is securely fastened by gently rotating it clockwise.

III. Product Operation:

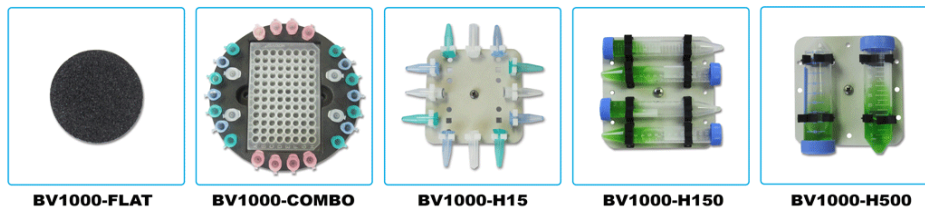
Set the mode switch to the desired setting of ON (continuous mixing) or TOUCH. In touch mode the tube to be mixed is pressed downward against the head to activate the vortex mixing action. Touch the tube to be mixed to the head and observe the vortex mixing action. Adjust the SPEED knob to obtain more or less vortex mixing action.

IV. Care and Maintenance:

No routine maintenance is required other than to keep the unit clean. Cleaning can be done with a damp cloth. Avoid the use of solvents as they may attack the rubber cup head or the ABS plastic housing.

V. Accessory Heads:

A wide variety of accessory heads are available including a Flat Head for multiple tube mixing, a Combo Head which holds up to 60 small tubes or one micro plate. Please contact your dealer if you need one of these accessory heads.



BV1000-FLAT

BV1000-COMBO

BV1000-H15

BV1000-H150

BV1000-H500

Website: www.BenchmarkScientific.com **E mail:** info@BenchmarkScientific.com

Tel: 908-769-5555

Fax: 908-222-1864

Rev. 8-11