



BenchRocker™ 2D Variable Speed/Angle Platform Rocker Model BR2000 Instruction Manual

Benchmark's BenchRocker™ 2D Platform Rocker is a quiet, high capacity, platform rocker whose speed and tilt angle are easily user adjustable. The BenchRocker™ 2D is a multi-purpose rocker serving most "see-saw" type rocking applications. It is ideal for blot washing and gel staining and de-staining. The BenchRocker™ has a 2.0kg (4.4lb) load capacity, comes with a flat, non-skid rubber mat. An optional stacking platform and dimpled mat are available accessories. The BenchRocker™ 2D will operate in cold rooms down to 4°C and in incubators up to 65°C.

I. Product Specifications:

Speed: User adjustable from 2rpm to 30rpm (4 tilts per minute to 60 tilts per minute)

Tilt angle: User adjustable from 0 to +/-30 degrees tilt

Motor: Electronically controlled brushless gear motor (no carbon brushes to replace)

Platform: 35.6 x 30.0 cm (14" x 12") with 3 tie down points on each side

Load Capacity: 2kg (4.4 lbs) centered and balanced on the platform

Dimensions: wxdxh 35.6x30.0x16.5 cm (14x12x6.5") Max tilt height is 24.1 cm (9.5")

Weight: 3.4kg (7.4lbs)

Mat: Flat, non-skid, black, autoclavable

Electrical: 115VAC, 60Hz, 0.24 amps (Model BR2000)

230VAC, 50Hz, 0.12 amps (Model BR2000-230) CE compliant

Operating environment: 4°C (cold room) to 65°C (incubator)

Warranty: 2 years

II. Product Set-Up:

Place the rocker on a clean, flat, stable surface. Make sure that the power switch is in the OFF (not illuminated) position. Set the unit to the desired tilt angle (see below). Plug the rocker into a properly grounded outlet. Check that nothing blocks the up and down rocking action of the platform. **Set the Speed Control knob to the lowest speed setting (full counter clock wise).**

III. Product Operation:

Load the platform from the center out, making sure the items placed on the platform create a balanced load around the center line of the platform. This is important as large, overhung platform loads may cause the rocker to tip over. Turn the unit on by activating the ON/OFF switch so that it is illuminated. Rotate the Speed Control knob clockwise to set the desired rocking speed.

IV. Tilt Angle Adjustment:

The tilt angle is adjusted at the back of the unit. First loosen the Safety Shield knob and move the Safety Shield aside to gain access to the tilt adjustment screw. Loosen the tilt adjustment screw with the 2.5mm hex Allen wrench provided with the unit. Sliding the black oval adjustment bar will increase or decrease the platform tilt angle. Once the platform angle is selected, re-tighten the tilt angle adjustment screw with the Allen wrench. **VERY IMPORTANT!** Position the safety shield so that it is centered over the black oval adjustment bar, then hand tighten the Safety Shield knob. The safety shield is used to keep fingers out of the rotating mechanism.

V. 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 damage the product housing.

website: www.BenchmarkScientific.com e mail: info@BenchmarkScientific.com