



BenchRocker™ 3D Variable Speed/Angle Platform Rocker/Rotator Model B3D2300 Instruction Manual

Benchmark's BenchRocker™ 3D Rocker Rotator is a quiet, high capacity, gyratory 3D rocker who's speed and tilt angle are easily user adjustable. The BenchRocker™ 3D produces a simultaneous rocking and rotating low speed action that is ideal for blot washing, gel staining and de-staining applications. At higher speeds, the BenchRocker™ 3D produces a vigorous gyratory action that is excellent for mixing sample tubes of all sizes. The BenchRocker™ 3D has a 2.0kg (4.4lb) load capacity and comes with both flat and dimpled rubber mats. An optional stacking platform is also available. The BenchRocker™ 3D 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

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

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

Platform: 300 x 300mm (12" x 12") with multiple tie down points on each side

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

Dimensions: wxdxh: 30.0 x 30.0 x 20.0 cm (12 x 12 x 8") Max tilt height is 29.2 cm (11.5")

Weight: 3.3kg (7.2lbs)

Mats: Both flat and dimpled mats are supplied and are autoclavable

Electrical: 115VAC, 60Hz, 0.24 amps (Model B3D2300)

230VAC, 50Hz, 0.12 amps (Model B3D2300-230) **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 rotating and up/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 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. Rotate the Speed Control knob clockwise to set the desired speed.

IV. Tilt Angle Adjustment:

The tilt angle is adjusted by hand. First remove the mat(s) from the platform. Then observe the hole at the center of the platform and the RED mark on the screw seen through the hole. This RED mark is the key to adjusting the platform tilt angle. **To increase the tilt angle**, grasp the platform and push down on the platform edge or corner pointed to by the red mark while simultaneously lifting up on the opposite platform edge or corner. The platform will loosen from its position allowing the tilt angle to be set by hand. Once the new tilt angle is selected simply replace the mat(s) on the platform. **To decrease the tilt angle**, grasp the platform and lift up on the edge or corner pointed to by the red mark while simultaneously pushing down on the opposite platform edge or corner.

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.