



BlotBoy™ and Mini BlotBoy™ Rockers Models B3D1308 and B3D1008 Instruction Manual

Benchmark's BlotBoy™ and Mini BlotBoy™ are three dimensional rocking mixers™ with their angle and speed optimally set for blot washing and gel staining / de-staining applications. The BlotBoy™ comes with a 30.0 x 30.0 cm (12 x 12") platform and the Mini BlotBoy™ comes with a 27.1 x 19.6 cm (10.5 x 7.5") platform. Both have flat, non-skid, rubber mats. The BlotBoy™ and Mini BlotBoy™ also make excellent general purpose lab rocker-rotators. Both have full 1.6kg (3.5 lb) load capacity. The BlotBoy™ and Mini BlotBoy™ 3D Rotators will operate in cold rooms down to 4°C and in incubators up to 65°C.

I. Product Specifications:

Speed: fixed at 12rpm (10 rpm for the 230 volt models)

Tilt angle: +/- 8 degrees from the horizontal

Motor: Long life AC gear motor (no DC motor carbon brushes to replace)

Platforms: BlotBoy™ 30.0x 30.0 cm (12 x 12") with 3 tie down points on each side

Mini BlotBoy™ 26.7 x 19.1 cm (10.5 x 7.5") with 3 tie down points each side

Mat: Flat, non-skid, black, autoclavable

Load Capacity: 1.6kg (3.5 lb) centered and balanced on the platform

Dimensions: BlotBoy™ (wxdxh) 30.0x30.0x21.1 cm (12x12x8.3") **Weight:** 2.0kg(4.4lbs)

Mini BlotBoy™ (wxdxh) 26.7x19.1x20.3 cm (10.5x7.5x8.0") **Weight:** 1.7kg(3.8lbs)

Electrical: 115VAC, 60Hz, 0.4 amps (Models B3D1308 and B3D1008)

230VAC, 50Hz, 0.2 amps (Models B3D1308-230 and B3D1008-230) **CE** compliant

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

Warranty: 2 years

II. Product Set-Up:

Place the rotator on a clean, flat, stable surface. Make sure that the power switch is in the OFF (not illuminated) position. Plug the rotator into a properly grounded outlet. Check that nothing blocks the up-down and rotating action of the platform.

III. Product Operation:

Load the platform from the center out, making sure the items placed on the platform form a balanced load around the center of the platform. This is important as large, overhung loads may cause the rotator to stall or tip. Turn the unit on by activating the ON/OFF switch so that it is illuminated. The rotator will begin to rotate.

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 mat or ABS plastic housing.

V. Accessories: Stacking Platforms and Mats:

Stacking platforms are available that allow doubling of the work area. There is a 7.0 cm (2.75") separation between the two platforms, allowing access to the content of both trays. The stacking platform comes with a flat mat. Additional dimpled and flat mats are also available.

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

Tel: 908-769-5555

Fax: 908-222-1864

Rev. 8-10