

Assurance® GDS Vortex Mixer

Assembly

For use with Assurance® GDS assays, remove the standard attachment (gently pull up at the corners) and snap the microtiter plate attachment onto the vortex mixer (align with top plate and press on side tabs of attachment).

Operation

- Place the GDS sample wells base on the vortex mixer's plate mounting attachment. Ensure that it is firmly secured to the mounting attachment. Ensure that the wells are covered with adhesive film before turning the shaker on.
- Turn the mixer on by pressing the POWER switch. To start, press MODE (to run).
- Turn the speed control knob slowly to set speed. See kit directions for speed settings.
- To stop the mixer, press MODE (to touch).

Important Safeguards

When using electrical equipment, basic safety precautions should always be followed to reduce the risk of fire, electric shock, and personal injury, including the following:

- Read all instructions before use.
- Do not immerse the mixer in water or liquids.
- To avoid electric shock, do not open the housing. Refer servicing to qualified personnel only.
- Be sure the power cord supplied with your mixer matches your AC power requirements.
- Do not operate if the equipment malfunctions or has been damaged in any manner.
- Use equipment only as specified in these directions.

Note: This product is for laboratory use only.

Cleaning

Unplug the mixer.

If contamination is suspected, moisten paper towel with bleach solution and wipe mixer with 10% bleach solution. Avoid spraying bleach solution directly onto surface. Allow bleach solution to remain on surfaces for a minimum of 15 min before wiping clean with 70% isopropyl alcohol solution.

To prepare 10% bleach solution add 10 mL of commercially available bleach containing at least 5% sodium hypochlorite to 90 mL of deionized water. The minimum final concentration of sodium hypochlorite in the bleach solution should be 0.5%. The bleach solution is stable for 7 days from preparation. To prepare 70% isopropyl alcohol solution add 70 mL of pure isopropyl alcohol to 30 mL of deionized water or buy commercial 70% isopropyl alcohol.

Caution: Follow the above listed safeguards when cleaning the instrument.

Manufacturing Entity

BioControl Systems, Inc, 12822 SE 32nd St, Bellevue, WA 98005, USA.

BioControl Systems, Inc is an affiliate of Merck KGaA, Darmstadt, Germany

Merck KGaA
Frankfurter Strasse 250
64293 Darmstadt
Germany
sigmaaldrich.com

