

Thermo Scientific

2000 mL SteriBIOS Bottle

Instructions for Use



Preface

Before starting to use the bottles read through this manual carefully and follow the instructions. Failure to follow the instructions and safety information in this manual will result in the expiration of the seller's warranty.

Intended Use

The 2000 mL SteriBIOS Bottle is intended for centrifugation in biotechnological applications. The bottle is specifically designed for use in combination with Sorvall BIOS 16 centrifuges and 8 x 2000 mL or 6 x 2000 mL swinging bucket rotors, respectively.

The 2000 mL SteriBIOS bottle is not approved or intended for, and should not be used for medical, clinical, surgical or other patient oriented applications.

Rotor Compartment Load

Bottles may be filled with up to 2000 mL. At 2000 mL the liquid level will be at the first thread of the bottle neck. Minimum fill level is 1000 mL. Refer to the drawing to see minimum fill level.

CAUTION If the maximum compartment mass of 2300 gram is exceeded, the speed must be reduced. Refer to centrifuge manual for reduction method.

Failure to do so can result in personal injury and/or centrifuge damage.

Chemical Resistance

The compatibility between chemicals and plastic centrifuge ware is affected by temperature, chemical concentration, g-force, length of run and other factors. Check the centrifuge manuals for chemical compatibility ratings. Because of the stresses associated with centrifugation, these ratings are a general guide only. We recommend a trial run before using a certain chemical.

Instruction

- CAUTION** Inspect the article upon delivery. Make sure that the protective PE bags are not damaged. If the PE bags are damaged, do not use the bottles, sterilization may be compromised.
- 2000 mL SteriBIOS Bottle requires its closure.
- Fill the bottle and place the closure onto the bottle, turn cap clockwise until hand-tight.
 - Before placing bottles into the rotor, weigh bottles to make sure that the bottles are balanced within the acceptable limit for the respective rotor being used. Consult the centrifuge manual for the acceptable imbalance tolerance.
 - After the run turn the closure counterclockwise to open the bottle.
- NOTICE** Before each use, inspect bottles for signs of wear including cracks, crazing, discoloration, yellowing, brittleness, deformation, surface abrasions or chemical attack. Immediately discard bottles showing any signs of wear.

Disposal

Do not recycle single use material. Follow your local regulations for disposal of waste and use proper protective equipment.

Product Specification

| Cat. No. | 75003871 |
|-------------------------------|---|
| Product Name | Thermo Scientific 2000 mL SteriBIOS Bottle |
| Description | 2000 mL bottle with closure; single-use, sterilized |
| Package | Set of 2 |
| Material Bottle | Polypropylene |
| Material Closure | Polypropylene |
| Max. Capacity [mL] | 2000 |
| Maximum Sample Density [g/mL] | 1.2 |
| Bottle Life Time | Single Use |
| Minimum Fill Level [mL] | 1000 |
| Temperature | In temperature to 4-22°C |
| G-force | In g-force to max. 7187 x g |
| Run Duration | If fill volume is between 1900-2000 mL, then no limitation If fill volume is between 1000-1899 mL, then limitation is max. 2 hours run duration |
| Centrifuge | Thermo Scientific Sorvall BIOS 16 |
| Rotor | 8 x 2000 mL 6 x 2000 mL |
| Autoclavable | N/A |
| Certifications | <ul style="list-style-type: none">Endotoxin per USP <85>Non-cytotoxic per USP <87>USP Class VI per USP <88>Particle Count Test per USP <788>BSE and TSE transmission safe |
| Sterilization | Sterilized using Beta irradiation following standards ISO 11137-1 and ISO 11137-2 |



Drawing

