

Guidelines for Use of BD Falcon™ Erlenmeyer Flasks

INTENDED USE

BD Falcon Erlenmeyer Flasks are disposable, reusable culture vessels that will allow cell or bacteria growth with uninhibited performance. Whether it is a cell biology, micro or molecular biology application BD Falcon Erlenmeyer Flasks are a proven, reliable, high quality tool that provides the flexibility of reuse or single use for when biohazardous or radioactive materials are used. The 2-in-1 cap is a patented feature allowing for vented or non-vented application with the same cap.

Erlenmeyer flasks made from glass have been used in the laboratory for over a century. As such, the shape and function of these containers has been standardized. They are ideal vessels for suspension culture and reagent storage. A polycarbonate (PC) flask that is provided sterile to an SAL 10^{-6} and can be autoclaved is an easier to use, safer product that is also certified pyrogen free and nuclease free.

STORAGE AND STABILITY

Store at room temperature (15-30° C).

PRODUCT SPECIFICATIONS

- Made from optically clear PC resin provides uninhibited visualization of sample or media
- 2-in-1 cap allows for vented or non-vented applications using the same product providing flexibility in their use
- Autoclavable(see instructions on reverse side) can be reused repeatedly; Media can be made up and autoclaved in same vessel
- Individually packaged, sterile, easy to use, maintains sample integrity
- Sterility to SAL 10^{-6} provides confidence in assay results especially during long term culture work
- Nuclease-Free for molecular biology or genomic work
- Non pyrogenic for cell culture
- Meets USP Class VI

RECOMMENDED CONDITIONS

Fill volume

The recommended fill volume is 25% to 40% of the nominal flask volume.

Seeding density

Start with a seeding density of 1×10^5 cells/ml. The optimal seeding density for your cell type with BD Falcon Erlenmeyer flasks will need to be optimized in order to achieve the best results. Results will vary depending on the cell type and media conditions.

Shaker speed

It is recommended to set the rotational speed between 50 rpm and 200 rpm. The optimal speed is cell type dependent. Optimization of the rotational speed should be performed to achieve the best results

STEAM STERILIZATION INSTRUCTIONS – see reverse side



STEAM STERILIZATION INSTRUCTIONS

The BD Falcon Erlenmeyer flasks are manufactured with polycarbonate that is designed to withstand repeated steam sterilization cycles. The following instructions are provided to aid in successfully using your labware as an ideal alternative to glass

For Filled Bottles or Flasks: Loosen the cap so it is barely engaged and sterilize according to the lab protocol. Allow the vessels and content to cool to 55°C or lower before securing the cap. Tightening caps on hot bottles will lead to distortion of the bottle as the contents cool down. The recommended sterilization temperature is 121°C at 15 psig, up to 60 minutes.

For Empty Containers: The caps should be removed from the vessels and sterilized in autoclave bags. As with glassware, cover the necks and tops of the flasks/bottles with aluminum foil. The foil should be crimped tightly enough so it does not fall off during autoclaving. A small piece of autoclave indicator tape should be used to secure the foil to the vessel, and to indicate the vessel has been exposed to steam.

Caution: The flasks may “implode” due to rapid exhaust setting (sudden change in temperature and pressure). For best results, use a slow exhaust cycle on the sterilizer.

DO NOT OVER STERILIZE. PC SHOULD NOT BE STERILIZED AT TEMPERATURES HIGHER THAN 129°C.

CUSTOMER AND TECHNICAL SERVICE

For technical assistance, contact Technical Service at:

Tel: 877.232.8995 or 978.901.7389 Fax: 978.901.7491; e-mail: Labware@bd.com

To place an order in the U.S., contact Customer Service at:

Tel: 877.232.8995 Fax: 800.325.9637 or 858.812.8889

Outside the U.S., contact your local distributor or nearest BD Biosciences office.

Visit our website www.bdbiosciences.com/discovery_labware for additional information

BD, BD logo, and all other trademarks are the property of Becton, Dickinson and Company. ©2006 BD