

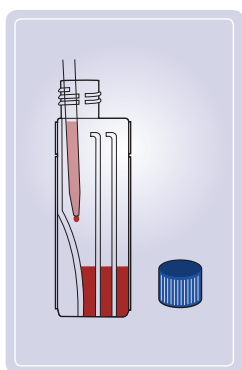
BD Falcon™ Cell Culture Multi-Flask

Guidelines for Use

INTRODUCTION

BD Falcon™ Cell Culture Multi-Flasks are available in 3- and 5-layer formats and provide 525 cm² or 875 cm² cell growth surface areas, respectively, which is equivalent to 3 or 5 times the surface area of a T-175 flask. BD Falcon Multi-Flasks offer the same benefits as the current format but in a more efficient and higher capacity design enabling you to grow more cells faster, easier, and in a more convenient manner. The footprint is the same as the T-175, so adaptation to your current equipment should be simple. This device allows you the benefit of preparing your cell suspension within the flask and removing your media or reagents using a pipet. This reduces the risk of contamination and speeds up the procedure.

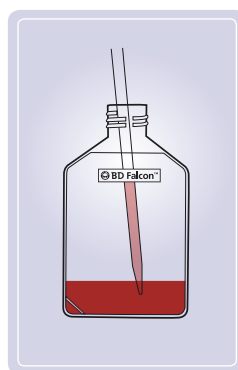
ADDING MEDIA AND PREPARING CELL SUSPENSION WITHIN THE BD FALCON MULTI-FLASK



1. Add required amount of medium into BD Falcon Multi-Flask by pipet or by pouring using typical culture volumes of 25-50 ml per layer.
2. To avoid foaming of medium, allow liquid stream to flow along the slope of the BD Logo side of the BD Falcon Multi-Flask.

Helpful hint: A 10 ml pipet allows media to be dispensed at the bottom of the vessel.

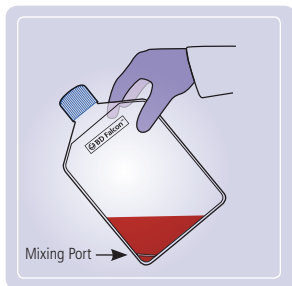
25-100 ml pipets allow media to be dispensed just past the BD Logo.



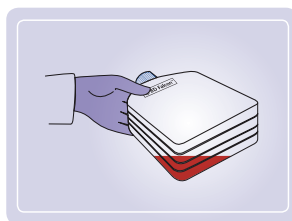
3. Dispense cell suspension from a concentrated stock into the growth medium using a BD Falcon 10 ml pipet (cat. no. 357551 or 356551). Be sure to dip the pipet tip into the medium.

Note: The seeding density will vary depending on the cell type, medium, and culture duration needs. Begin with the same seeding density on a cells per centimeter square area to that used in standard flask for the cell type used.

MIXING OF CELLS IN BD FALCON MULTI-FLASK



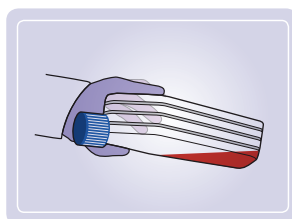
4. Mix position: Hold the BD Falcon Multi-Flask upright with the BD Logo facing you and tilt counter-clockwise to a 45° angle.



- 5a. Holding at the same angle, gently rotate the BD Falcon Multi-Flask forward (neck pointing away from you).

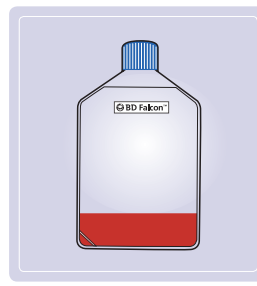
- 5b. Then, gently rotate it backward (neck pointing towards you).

Note: With each tilt, hold until liquid in the top layer drains fully.



6. Repeat Step 5 to ensure proper mixing. Bring back to mix position, as shown in Step 4. Then, proceed to Step 7 for equilibration.

EQUILIBRATING FLUID



7. After mixing or adding cell suspension, place the BD Falcon Multi-Flask vertically on a flat work surface to equalize liquid volume among all the layers.

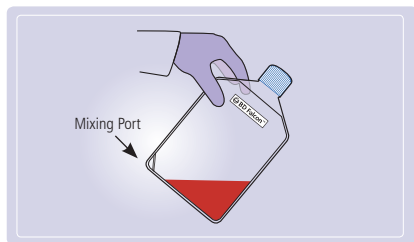
ALTERNATE PROTOCOL

Adding cell suspension prepared external to the BD Falcon Multi-Flask

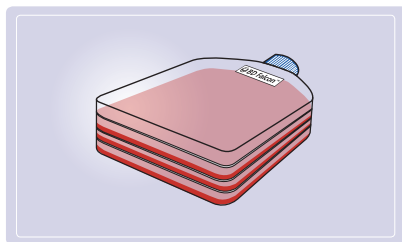
1. Create cell suspension externally from the BD Falcon Multi-Flask.
2. Add required amount of cell suspension into the BD Falcon Multi-Flask by pipet or by pouring.
3. To avoid foaming of medium, allow liquid stream to flow along the slope of the BD Logo side of the BD Falcon Multi-Flask.



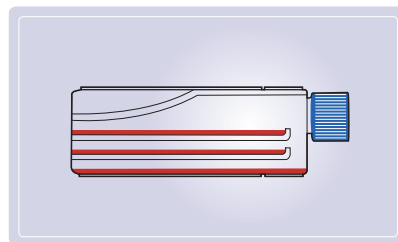
PARTITION AND DISTRIBUTE LIQUID ONTO EACH LAYER



8. Hold the BD Falcon™ Multi-Flask upright with the BD Logo facing you and tilt clockwise to a 45° angle on a flat work surface to partition the liquid into each of the layers. This position is recommended for transport.



9. While holding the BD Falcon Multi-Flask at a 45° angle, gently lay it flat onto the work surface with BD Logo facing up.

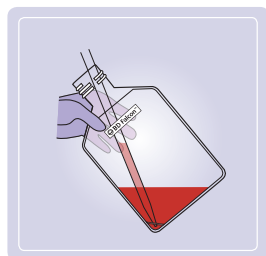


10. After placing the BD Falcon Multi-Flask flat on a work surface, gently rock back and forth and side-to-side to distribute cells evenly onto culture surfaces – taking care not to spill liquid from each layer.

MEDIA REMOVAL

If exchange of media is required, follow Steps 7-10.
You may choose to either aspirate or pour the media from the BD Falcon Multi-Flask.

ASPIRATING METHOD

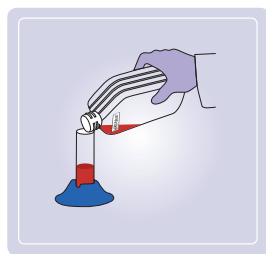


11. To aspirate or remove media tilt the BD Falcon Multi-Flask, with the BD Logo facing you, counter-clockwise to a 45° angle while inverting the multi-flask toward you.

12. Then, tilt the BD Falcon Multi-Flask to the right, continuing to aspirate all residual media.

Helpful hint: Aspirate media using a BD Falcon 2 ml aspirating pipet (cat. no. 357558).

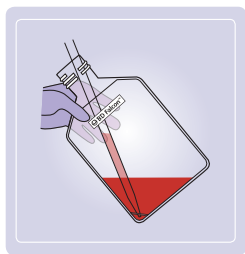
POURING METHOD



13. With BD Logo facing down, pour spent media from the BD Falcon Multi-Flask.

Helpful hint: Pouring is easier when the BD Logo is facing you and the multi-flask is tilted at a counter-clockwise 45° angle.

CELL HARVESTING



14. Add dissociating reagent (≥ 5 ml per layer) based on preferred protocol and bring to mix position (Step 4). Then, follow Steps 7-10.

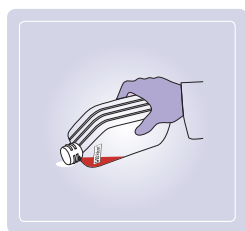
15. Neutralize with inactivating solution and mix following Steps 4-10. Gently swirl to dislodge cells completely.

16. Pipetting Method: Follow “Media Removal” protocol but collect cell suspension using a BD Falcon 10 ml serological pipet (cat. no. 357551).

17. Follow Step 13 “Pouring Method”. Pour detached cell suspension into a BD Falcon conical tube (cat. no. 352070).

18. Rinse with additional media as needed.

Helpful hint: Pouring is easier when the BD Logo is facing you and the multi-flask is tilted at a counter-clockwise 45° angle.



PRODUCT SPECIFICATIONS

WORKING VOLUME RANGE	≥ 5 ml per layer for dissociating ≤ 50 ml per layer for cell expansion
MOLDED-IN GRADUATIONS	0 to 50 ml per layer in 10 ml increments
GRADUATION ACCURACY	10%
CAP VENT MEMBRANE	0.2 μ m hydrophobic membrane
CELL GROWTH SURFACE	Tissue culture-treated, optically clear

ORDER INFORMATION

BD Falcon Cell Culture Multi-Flask

Description	Qty./Pack	Qty./Case	Cat. No.
3-LAYER TISSUE CULTURE-TREATED, 525 CM ²	2	12	353143
5-LAYER TISSUE CULTURE-TREATED, 875 CM ²	1	8	353144

U.S. Orders

tel: 877.232.8995
e-mail: BDBCcustomerService@bd.com

Purchase Orders should be made out to:
BD Biosciences
2350 Qume Drive
San Jose, CA 95131

Outside the U.S.

Contact your local distributor or visit bdbiosciences.com/offices to locate your nearest BD Biosciences office.

ACCESSORIES

Description	Qty./Pack	Qty./Case	Cat. No.
BD Falcon Pipets			
ASPIRATING PIPET 2 ML INDIVIDUALLY WRAPPED	50	200	357558
10 ML INDIVIDUALLY WRAPPED PAPER/PLASTIC	50	200	357551
10 ML INDIVIDUALLY WRAPPED ALL-PLASTIC	50	200	356551

BD Falcon Tubes

50 ML POLYPROPYLENE CONICAL TUBE	25	500	352070
175 ML POLYPROPYLENE CONICAL TUBE	8	48	352076
225 ML POLYPROPYLENE CONICAL TUBE	8	48	352075

U.S. Orders

Contact your authorized distributor to place your order.



BD Biosciences

Two Oak Park
Bedford, MA 01730
tel: 877.232.8995
fax: 800.325.9637
bdbiosciences.com