

BD Biosciences - Discovery Labware
296 Concord Road
Billerica, MA 01821
tel: 877.232.8995
fax: 800.325.9637
bdbiosciences.com



March 2007

Dear Valued Customer:

We are writing to notify you that **BD Biosciences** – Discovery Labware is implementing a packaging enhancement for our **BD Falcon™ 15 ml Conical Tubes**. Our records indicate you have purchased these products within the last two years. To better serve you, we have changed the packaging design; however, the primary packaging material will remain high-density polyethylene. The case count will remain 500 tubes per case; however, the shelf-pack count has been reduced from 125 to 50 tubes. We're confident you'll find the change will enable you to more easily stack, open, and use our tubes for all your lab needs.

This change impacts the following products:

- BD Falcon 15 ml Polystyrene Conical Tube (Cat. No. 352095)
- BD Falcon 15 ml High-clarity Polypropylene Conical Tube (Cat. No. 352096)
- BD Falcon 15 ml High-clarity Polypropylene Conical Tube w/Rack (Cat. No. 352196)

Measurement changes to the case dimensions are as follows:

Cat. No.	Current Case	New Case
352095	12.8" x 11.3" x 10.6"	13.2" x 11.07" x 11.12"
352096	12.8" x 11.3" x 10.6"	13.2" x 11.07" x 11.12"
352196	27.5" x 12.7" x 6.0"	13.2" x 11.07" x 13.12"

Please note that for the first two months of production with the new packaging, a product insert will accompany the product, detailing the changes.

If you would like more information about this change, you may contact your local BD Biosciences representative or BD Biosciences Technical Support at **877.232.8995**. Outside the U.S., visit our Website at **bdbiosciences.com/offices** to locate an office near you.

We remain committed to serving all your research needs, and appreciate your business. If we may better serve you in any way, please contact us.

Sincerely,

Heather Donovan
Quality Systems & Compliance Manager
heather_donovan@bd.com

Patricia Daggett
Product Manager, Fluid Handling and TC Products
patricia_daggett@bd.com