

**BD Biosciences**  
Discovery Labware  
296 Concord Road  
Billerica, MA 01821 USA  
tel: 978.671.1388  
fax: 978.901.7490  
bdbiosciences.com



September 2007

Dear Valued Customer:

**BD Biosciences** – Discovery Labware is implementing a manufacturing location and a case dimension change for our BD Falcon™ 60 mm Easy-Grip Cell Culture Dish, **Cat. Nos. 353004** (Tissue Culture-Treated) and **351016** (Non Tissue Culture-Treated). The manufacturing location has changed from the Plymouth, U.K. facility to the Durham, N.C. facility. Please note that both of these BD Biosciences facilities manufacture under the ISO 9001:2000 Standard, and the current FDA Quality System Regulation in Chapter 1, Title 21, Section 820 of the Code of USA Federal Regulations.

**Measurement changes to the case dimensions:**

Current case	New case
12 3/8" x 9 7/8" x 15 3/8"	12 3/8" x 10" x 15 1/4"

For the first 3 months of sales, product manufactured in the Durham, N.C. location will be identified by a case insert. They will also be permanently identifiable by the Durham, N.C. address shown on the packaging.

If you would like more information about this change, please contact your local BD Biosciences representative or BD Biosciences Technical Support at **877.232.8995** or [labware@bd.com](mailto:labware@bd.com) (U.S.). For assistance outside the U.S., please reference your BD Biosciences catalog for the phone number of the BD office nearest you or visit our website at [bdbiosciences.com](http://bdbiosciences.com), click "Support", then "Contact BD" for up-to-date worldwide contact information.

Thank you for your loyal business with BD Biosciences.

Sincerely,

A handwritten signature in black ink, appearing to read 'Heather Donovan', with a long horizontal flourish extending to the right.

Heather Donovan  
Quality Systems & Compliance Manager

A handwritten signature in black ink, appearing to read 'William R. Beltz', with a long horizontal flourish extending to the right.

William R. (Bob) Beltz, Ph.D.  
Product Manager