

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 the following modifications to our BD Falcon™ Petri Dish 100 mm x 15 mm, Cat. No. 351029:

— **New Resin Supplier** - The resin used in the manufacture of the dish remains crystal-grade polystyrene however we have validated an alternative supplier.

— **Case Measurement** -

Current case	New case
19.3" x 19.3" x 12.9"	21.5" x 20.2" x 14"

Note: The case count will remain 500 dishes per case and 20 dishes per sterile package.

— **Product Dimension Change** - Less than 1% dimensional changes to the dish and lid.

Materials manufactured under this change notification will be identified by a sticker on the case stating "**New and Improved Packaging**", for a minimum of the first 3 production lots.

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 (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, click "Support", then "Contact BD" for up-to-date worldwide contact information.

Thank you for your loyal business with BD Biosciences.

Sincerely,

Heather Donovan
Quality Systems & Compliance Manager

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