

## New Y-Fitting for the Model 475 Gradient Delivery System

The Y-fitting is used with the Gradient Delivery System to mix the high and low density gel solutions during casting of gradient gels. This new Y-fitting is a replacement for the existing Y-fitting. There are no metal tubing or screws associated with the new Y-fitting. The new Y-fitting will generate the same linear gradients as the old Y-fitting.

**DCode™ Universal Mutation Detection System customers do not need to make any changes to their DCode instruction manual.**

Please make the following new changes to your *D GENE* and Model 475 Gradient Delivery System instruction manuals.

### Changes for the *D GENE* Instruction Manual

Section 6.3, step 3, page 19, Disregard the first sentence and replace the second sentence with "Connect the free end of the 9 cm Tygon tubing to one end of the Y-fitting and connect a luer coupling to the other end of the Tygon tubing."

Section 6.3, step 10, page 20, Replace the first sentence with "Attach the tubing from the high density syringe to one end of the Y-fitting."

Section 6.3, step 14, page 20, Disregard the third sentence "Detach the syringe tubing from the metal tubing on the Y-fitting."

Section 6.4, step 3, page 21, Replace the second sentence with "Connect the free end of the 9 cm Tygon tubing to one end of the Y-fitting and connect a luer coupling to the other end of the Tygon tubing."

Section 6.4, step 6, page 22, Replace the first sentence with "Attach the tubing from the high density syringe to one end of the Y-fitting."

Section 6.4, step 9, page 22, Disregard the third sentence "Detach the syringe tubing from the metal tubing on the Y-fitting."

### Changes for the Model 475 Gradient Delivery System Instruction Manual

Section 1.2.10, page 3, Disregard the second sentence "Three metal tube fittings are inserted into the Y-fitting and held in place by three screws."

Section 2.1, step 2, page 4, Replace the second sentence with "Connect the free end of the 9 cm Tygon tubing to one end of the Y-fitting and connect a luer coupling to the other end of the Tygon tubing."

Section 2.1, step 8, page 4, Replace the fifth sentence with "Attach the tubing from the high density syringe to one end of the Y-fitting."

Section 2.1, step 13, page 5, Disregard the third sentence "Detach the syringe tubings from the metal tube on the Y-fitting."

Section 2.2, step 2, page 6, Disregard the first sentence and replace the second sentence with "Connect the free end of the 9 cm Tygon tubing to one end of the Y-fitting and connect a luer coupling to the other end of the Tygon tubing."

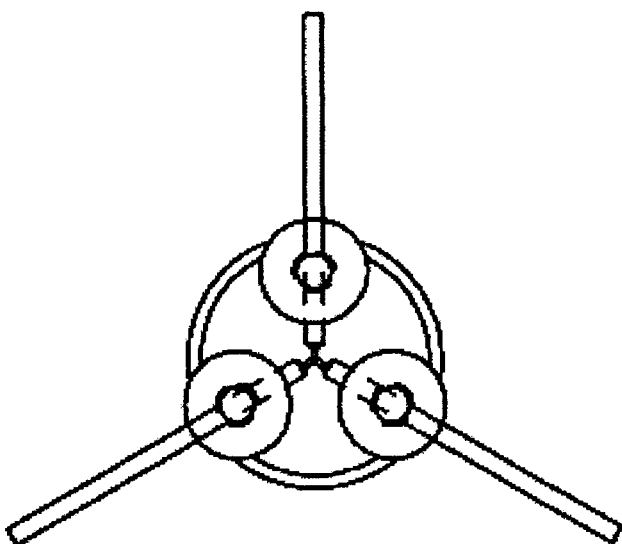
Section 2.2, step 4, page 6, Replace the fifth sentence with "Attach the tubing from the high density syringe to one end of the Y-fitting."

Section 2.2, step 8, page 7, Disregard the third sentence "Detach the syringe tubings from the metal tube on the Y-fitting."

Section 3, page 8, Disregard the third sentence "Detach the syringe tubings from the metal tubings on the Y-fitting."

We recommend these instructions be inserted into the instruction manual.

If you have any questions or would like further information, contact your local Bio-Rad representative, or in the US call 1-800-4BIORAD.



**Old Y-fitting**



**New Y-fitting**