



## Instruction for Converting a Phosphor Screen for Bio-Rad FX Systems

March 2001

1. Carefully remove the screen from the aluminum backing plate. Use a sharp knife or razor blade to lift a corner of the screen. Peel the screen up without bending it sharply, which can cause creases. If you have difficulty removing the screen, heat the glue with a hair dryer or place the whole screen into an oven set to warm (approximately 150 to 200 degrees F), and then try the above procedure.
2. Once the screen has been removed, place the screen face (white side) down onto a clean, dry surface (Caution: avoid water on the surface, since water can damage the screen permanently).
3. Peel off the adhesive backing on the screen backing about one third of the way and only cut off the paper backing on this portion of the screen. Place the exposed adhesive region of the screen so that it is centered on the screen from side to side and leaves a border of about 4-5 mm on the top of the screen (the screen backing is slightly smaller than the screen).
4. Smooth the screen backing with your hand to remove any air bubbles. Lift up the unattached portion of the screen backing and remove the rest of the adhesive backing from the screen backing. Smooth the screen backing onto the screen to remove any air bubbles as before.
5. Your screen is now ready to use in the Bio-Rad Molecular Imager FX systems.