



Instructions for Anode Plate Installation

Enclosed is your new anode plate electrode which has been redesigned to provide a more reliable electrical connection. Handle the plate at the edges or on the flange to maintain the clean appearance. If you have a recent model of the Semi-Dry blotter already equipped with the new electrode plate, simply replace the old plate making the same connections as on the original arrangement. If you own an older version of the unit, the anode electrode is a flat plate. You will notice the replacement electrode is different in appearance and will now extend over the edge of the spring-supported platform when assembled. Please follow the assembly instructions below to replace with the new electrode.

1. Remove the old electrode by unscrewing the four plastic screws which secure it to the platform.
2. Lift the old electrode plate off of the platform and discard.
3. Wipe off the foil surface. Steel wool may be used to clean off deposits if necessary. It is not necessary to remove the stainless steel foil although it will no longer be functional.
4. Unscrew the four white posts which hold the spring-supported platform in place.
5. Lift out the platform and detach the wire from the platform by unscrewing the metal screw.
6. Assemble the new electrode plate onto the platform and secure with the four plastic screws.
7. Anchor the wire from the base and the plate electrode together to the platform using the metal screw.
8. Place the assembled platform back onto the base making sure that the four springs remain upright.
9. Slide the four white posts back into the holes in the platform and screw them back into the base completing the assembly.

If you require further assistance, in the U.S. please telephone our Technical Service Department at 1-800-4BIORAD (1-800-424-6723). Outside the U.S., please contact your local Bio-Rad representative.