

**BIO-RAD**

### **Trans-Blot Plus Gel/Cassette Assembly Tray**

This tray is intended for use in blotting cassette assembly and gel staining. The lid imparts stability to the sides and lip of the tray. When liquid, gels or other materials are placed within the tray, it should not be transported without the lid in place.

*Bio-Rad Laboratories, 2000 Alfred Nobel Drive, Hercules, CA 94547*

4106243 Rev A

**BIO-RAD**

### **Trans-Blot Plus Gel/Cassette Assembly Tray**

This tray is intended for use in blotting cassette assembly and gel staining. The lid imparts stability to the sides and lip of the tray. When liquid, gels or other materials are placed within the tray, it should not be transported without the lid in place.

*Bio-Rad Laboratories, 2000 Alfred Nobel Drive, Hercules, CA 94547*

4106243 Rev A

**BIO-RAD**

### **Trans-Blot Plus Gel/Cassette Assembly Tray**

This tray is intended for use in blotting cassette assembly and gel staining. The lid imparts stability to the sides and lip of the tray. When liquid, gels or other materials are placed within the tray, it should not be transported without the lid in place.

*Bio-Rad Laboratories, 2000 Alfred Nobel Drive, Hercules, CA 94547*

4106243 Rev A

**BIO-RAD**

### **Trans-Blot Plus Gel/Cassette Assembly Tray**

This tray is intended for use in blotting cassette assembly and gel staining. The lid imparts stability to the sides and lip of the tray. When liquid, gels or other materials are placed within the tray, it should not be transported without the lid in place.

*Bio-Rad Laboratories, 2000 Alfred Nobel Drive, Hercules, CA 94547*

4106243 Rev A