



## *Important Notice*

### **Trans-Blot<sup>®</sup> Super Cooling Coil Catalog 170-3912**

1. Before using, check coil for any shipping/handling damage such as scrapes or abrasions. If damage is found, contact Bio-Rad Laboratories to arrange for replacement. Do not use a damaged coil.
2. This coil has been specifically designed for use with the Trans-Blot cell. Do not use in applications over 500 V DC.

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