
Installation Qualification and Operational Qualification Protocols for the PowerPac Universal Power Supply

**Catalog Number
164-5068**

(Part Number 4506017, Rev A)

BIO-RAD