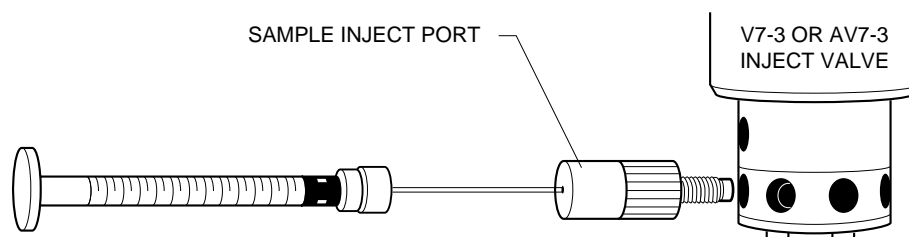


## **SAMPLE INJECT PORT**

**CATALOG NUMBER 750-0471**

### **To install the Sample Inject Port:**

1. Attach a 1ml disposable syringe to the 22 gauge (0.028" diameter) blunt tip injection needle supplied with the fittings kit. Place the ferrule end of the Injection Port against a clean flat surface and firmly push the needle through to the flat surface. The tip of the needle should be visible from the ferrule end. This initial fit is tight but will loosen with use.
2. With the needle still in the Sample Inject Port, screw the entire assembly into port 5 of the inject valve (V7-3 or AV7-3). Finger tighten until snug. Do not use a wrench.
3. Remove the syringe and needle.



*Installing the Sample Inject Port*

### **Technical Assistance**

For additional help, contact your local Bio-Rad representative. In the United States, call Technical Service at 1-800-4BIORAD.

### **Ordering Information**

750-0471 Sample Inject Port (1)