

## **Model 2110 Fraction Collector Sole Source Specifications**

The Model 2110 Fraction Collector includes the fraction collector base unit, carousel tube holder, 2-way stopcock, and user manual.

### **The Model 2110 Fraction Collector is capable of:**

- Time or drop collection modes (or volume collection mode, when connected to Model EP-1 Econo pump, BioLogic LP, or BioLogic DuoFlow system).
- Time intervals of .05–99.9 minutes.
- Small chromatography columns or gravity columns mount directly to the drop-forming arm to minimize dead volume.
- Silicone drop former accommodates either 1/8" or 1/16" tubing without the need for luer or ¼-28 fittings.
- Manually adjustable drop arm detector signal level.
- Optional Instrument Dust cover intended for use when the drop arm is in collect position. The inlet tube may be directed out the top opening or through the small opening at the bottom/rear of the instrument.
- Programmable and controlled by the BioLogic LP System. When used with the BioLogic LP system up to 50 collection windows are permitted per stored method.
- Controlled via the DuoFlow chromatography system software.
- Fraction event marks can be sent to either a chart recorder and recorded or event marks can be sent through the BioLogic LP controller to be recorded on the LP Data View collection software.
- The Model 2110 fraction collector meets I.E.C. 1010 safety certification, an international safety standard, and C.S.A. 22.2.