

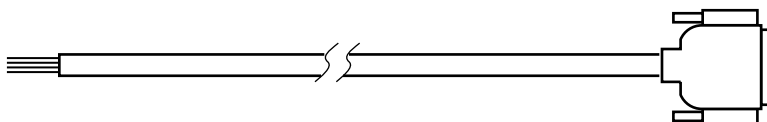
SYSTEM CABLE 5 FOR THE MODEL 2110 FRACTION COLLECTOR

CATALOG NUMBER 731-8265

The Model 2110 fraction collector can be connected to the BioLogic System using System Cable 5, a breakout cable which connects to the DB-9 connector on the Model 2110 Fraction Collector and to the AUX connector on the BioLogic Workstation. To connect to the BioLogic AUX connector, make the following connections:

- Black wire to AUX connector #5
- White wire to AUX connector #9

Cut all other wires as close to the shielding as possible and cover the ends with electrical tape



DB-9 Connector Pin number	Wire Color	Model 2110 Function	Signal Type
Pin 1	Black wire	External advance	Input TTL (active low)
Pin 2	Orange wire	Drops	Output TTL (active low)
Pin 3	Yellow wire	Advanced	Output TTL (active low)
Pin 4	Red wire	Ready	Output TTL (active low)
Pin 5	Brown wire	Alarm	Output TTL (active low)
Pin 6	Purple wire	External start	Input TTL (active low)
Pin 7	Blue wire	Advanced Contact 1	
Pin 8	Green wire	Advanced Contact 2	
Pin 9	White wire	Logic Ground	

Technical Assistance

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