

# Instructions for Proper Assembly of the PROTEAN<sup>®</sup> plus Dodeca<sup>™</sup> Cell Buffer Recirculation Path

1. Remove the PROTEAN plus Dodeca Cell from the shipping box and place it on the lab bench. Position the cell so that you can easily access the plumbing on the back of the instrument.
2. Remove the buffer recirculation pump from the shipping box. Place it on the same lab bench alongside the PROTEAN plus Dodeca Cell, positioned so that the front of the pump is facing you.
3. Install the Buffer Exhaust Tube to the lid of the cell. The connector is a quick-connect fitting. Push it into the lid connector until an audible click is heard.
4. Connect the outlet of the Buffer Exhaust Tube to the inlet of the Buffer Recirculation Pump.
5. Connect the outlet of the Buffer Recirculation Pump to the Buffer Recirculation Pump path at the bottom of the cell. The connector is a quick - connect fitting.

**NOTE THE BUFFER PATH. BUFFER IS PULLED FROM THE TOP OF THE CELL AND REINTRODUCED TO THE BOTTOM OF THE CELL.**

