



Rotofor® Preparative IEF Cell Heat Exchanger O-ring Kit

Catalog Number 170-2954

Replacing Heat Exchanger O-rings

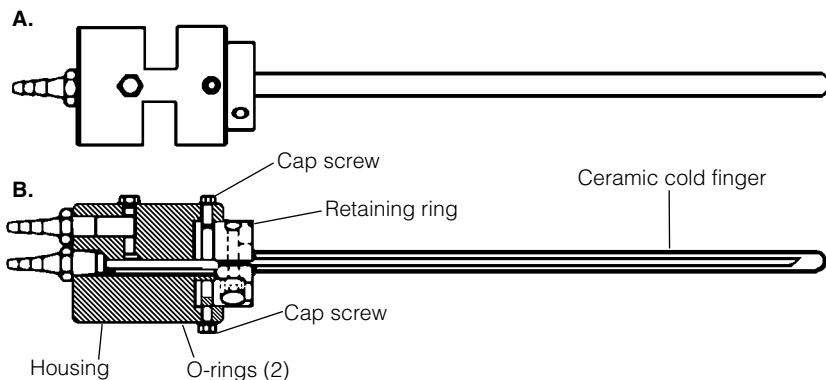


Fig. 1. A) Heat exchanger assembly. B) Section view of the heat exchanger assembly.

1. Remove the heat exchanger assembly from the Rotofor cell by disconnecting the coolant tubing. Remove the focusing and buffer chambers and the membrane core.
2. Loosen the cap screws by rotating them six complete turns counterclockwise. See Figure 1.
3. Remove the ceramic cold finger with retaining ring from the housing. Do not separate the ceramic cold finger from the retaining ring.
4. Remove the two worn O-rings. Wipe excess lubricant from housing and ceramic cold finger.
5. Install two new O-rings on ceramic cold finger and pack new O-rings with O-ring lubricant (supplied with the Heat Exchanger O-ring Kit).
6. Insert the ceramic cold finger with retaining ring and new O-rings in the housing.
7. Tighten the cap screws by rotating them six complete turns clockwise. Make sure the ceramic cold finger can rotate freely after the cap screws have been tightened.