

## CARBONATE REMOVAL DEVICE 300 (CRD-300) QUICKSTART

The CRD-300 must be handled with care to ensure proper operation. Fittings only need to be finger tightened.



**WARNING**

*Do not attempt to disassemble the CRD-300; it may result in irreversible damage.*

### 1.1. Hydrating the CRD-300

#### Step 1

Using a 5 cc disposable plastic syringe (P/N 016640) and the 10-32 Luer adaptor (P/N 046888), push approximately 3 mL of degassed DI water through the ELUENT IN port. Using a 5 cc disposable plastic syringe (P/N 016640) and the 1/4-28 Luer adaptor (P/N 024305), push 5 mL of degassed DI water through the REGEN IN port (Figure 1).

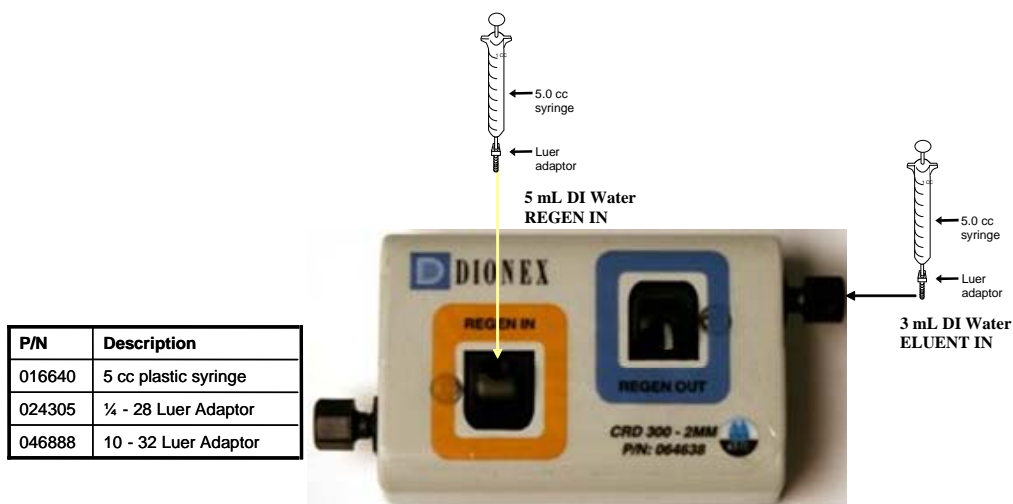


**NOTE**

*Step 1 can be accomplished by installing the CRD 300 in the system and connecting the ELUENT OUT port to the REGEN IN port on the CRD by using suitable tubing and pumping 5 mL of deionized water through the CRD ELUENT IN port. In the above step, it is recommended to bypass the guard and analytical columns.*

#### Step 2

Allow the CRD-300 to sit for approximately 10 minutes to fully hydrate the CRD-300 membrane.



**Figure 1**  
**Hydrating the CRD-300**

### 1.2. Backpressure Instructions



**WARNING**

*Total backpressure amounts exceeding >100 psi may cause irreversible damage to the CRD 300.*

The total backpressure to the suppressor eluent channel should be less than 60 psi. This includes the CRD 300, the cell, and the backpressure coil. Trim the backpressure coil if required to achieve <60 psi total backpressure. Refer to Section 3.4 of the CRD 300 Manual (P/N 065213) for backpressure measurement instructions.