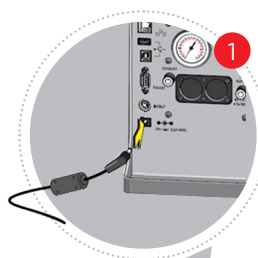


Switching Test Fluids

You must perform one of the following procedures when switching the fluids.


- Switching Test Fluids from Oil to Water
- Switching Test Fluids from Water to Oil

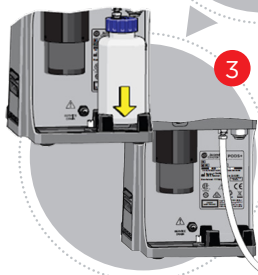
Switching Test Fluids from Oil to Water



1 Plug the external power supply jack into the PODS+ at the power port location.



2 Press the  key to power on the instrument.




3 Make sure the waste bottle is attached to the unit or drain tubing is connected to the drain port.

4 Fill the sample bottle with clean mineral spirits.









5 Place the sample bottle with clean mineral spirits in the Sample Cup Adapter and connect it to the instrument.

From the PODS+ Home screen, press the menu button next to  to display the Cleaning Routines Screen.



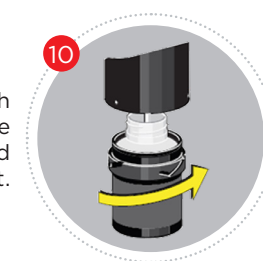
7 Setup the Cleaning Routine.







- Press the menu button next to .
- Press  on the keypad and select 50mL/min flow rate.
- Press  on the keypad and enter 55 for the volume.
- Press the menu button next to .
- Press the menu button to .

8 Press the menu button next to  to run the cleaning routine.

9 After sample run is complete, remove the sample bottle and re-fill the bottle with clean Isopropyl alcohol.








Place the sample bottle with clean Isopropyl alcohol in the sample cup adapter and connect it to the instrument.



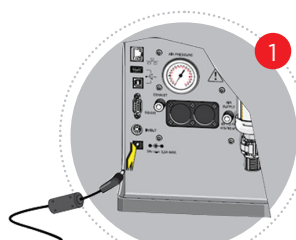
- 11 Setup the Cleaning Routine.
- A. Press the menu button next to .
 - B. Press  on the keypad and select 50mL/min flow rate.
 - C. Press  on the keypad and enter 55 for the volume.
 - D. Press the menu button next to .
 - E. Press the menu button to .
- 12 Press the menu button next to  to run the cleaning routine.
- 13 After sample is complete, remove the sample bottle and re-fill the bottle with clean DI water.



Place the sample bottle with clean DI water in the Sample Cup Adapter and connect it to the instrument.


- 15 Setup the Cleaning Routine.
- A. Press the menu button next to .
 - B. Press  on the keypad and select 50mL/min flow rate.
 - C. Press  on the keypad and enter 55 for the volume.
 - D. Press the menu button next to .
 - E. Press the menu button to .
- 16 Press the menu button next to  to run the cleaning routine.
- 17 When the sample run is complete, Press  to return to the Home screen.

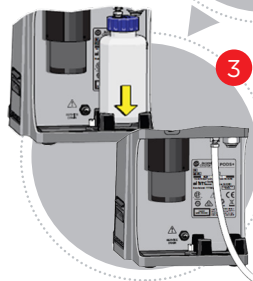
Switching Test Fluids from Water to Oil



1 Plug the external power supply jack into the PODS+ at the power port location.



2 Press the  key to power on the instrument.




3 Make sure the waste bottle is attached to the unit or drain tubing is connected to the drain port.

4 Fill the bottle with clean Isopropyl alcohol.









5 Place the sample bottle with clean Isopropyl alcohol in the Sample Cup Adapter and connect it to the instrument.



6 From the PODS+ Home screen, press the menu button next to  to display the Cleaning Routines Screen.

7 Setup the Cleaning Routine.

- A. Press the menu button next to .
- B. Press  on the keypad and select 50mL/min flow rate.
- C. Press  on the keypad and enter 55 for the volume.
- D. Press the menu button next to .
- E. Press the menu button to .






8 Press the menu button next to  to run the cleaning routine.


9 After sample is complete, remove the sample bottle and re-fill the bottle with clean Mineral Spirits.



Place the sample bottle with clean Mineral Spirits in the sample cup adapter and connect it to the instrument.

11 Setup the Cleaning Routine.









- A. Press the menu button next to .
- B. Press  on the keypad and select 50mL/min flow rate.
- C. Press  on the keypad and enter 55 for the volume.
- D. Press the menu button next to .
- E. Press the menu button to .

12 Press the menu button next to  to run the cleaning routine.

13 After sample is complete, remove the sample bottle and re-fill the bottle with clean DI water.



Place the sample bottle with clean MIL-5606 water in the Sample Cup Adapter and connect it to the instrument.

- 15 Setup the Cleaning Routine.
- A. Press the menu button next to .
 - B. Press  on the keypad and select 50mL/min flow rate.
 - C. Press  on the keypad and enter 55 for the volume.
 - D. Press the menu button next to .
 - E. Press the menu button to .
- 16 Press the menu button next to  to run the cleaning routine.
- 17 When the sample run is complete, Press  to run the cleaning routine one more time.
- 18 When the sample is complete, Press  to return to the Home screen.