

Best Practices

Solvents:

- Use clean borosilicate bottles only
- Prepare solvent volume to be used up within 1 – 2 days
- Use only HPLC-grade or better solvents filtered through 0.2 μm filters
- Label bottles correctly
- Reduce risk of exposure to direct sun light by using brown bottles

Powering up the pump:

- Use fresh solvents and solvent bottles
- Inspect solvent bottles and inlet filters for damages or coloring
- Check function and settings of the seal wash

Preparing the pump:

- Purge the pump to remove old solvent, salt deposits or particulate matter
- Condition the system with appropriate conditioning method

Daily tasks:

- Replace water based mobile phases daily
- Replace organic mobile phases every 2nd day
- Check seal wash solvent
- Use the Conditioning function with composition of your application

Weekly tasks:

- Change seal wash solvent and bottle
- Inspect solvent filters
- Flush all channels to remove salt deposits and particulate matter
- Check system backpressure and change filters if necessary

Pump shut-down:

- Flush all channels to remove salt deposits and particulate matter
- Flush the system with appropriate storage solvent
- Power down the system



Seal Wash (usage mandatory when installed)

Benefits of Seal Wash Operation:

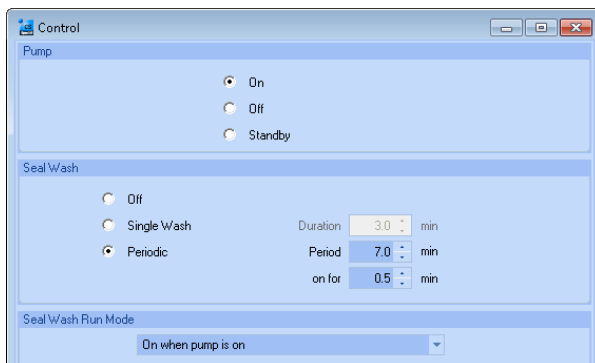
- Removal of particles and salt crystals, which might potentially damage the seals
- Lubrication of seal/piston interface
- Cooling of pistons

Seal Wash Operation:

- PERIODIC operation, e.g. 0.5 min every 7 min
- Use 10 % 2-Propanol in water (100 % 2-Propanol for normal phase applications)
- Position wash solvent bottle above and waste bottle below instrument
- Exchange solvent weekly. Do not recycle seal wash solvent
- Test Peristaltic Pump

Seal Wash Dialog in OpenLab Software (1290 Infinity Pumps only):

- Seal Wash settings are NO method parameters
- Find dialog under CONTROL
- On ERROR – seal wash is switched off
- At Power Off – seal wash is switched off



Seal Wash in 1290 Infinity II Pumps:

- PERIODIC operation for 0.5 min every 7 min when pump is switched on
- Controlled by the seal wash sensor
- Guided out to the waste on the right side



01200-90091

Part Number:
01200-90091

Edition: 04/2015
Printed in Germany

© Agilent Technologies, Inc 2015

Agilent Technologies, Inc
Hewlett-Packard-Strasse 8
76337 Waldbronn, Germany