

# 1290 Infinity II Bio and 1260 Infinity II Prime Bio LC

## Quick Reference Sheet

### Best Practices

Solvents:

- Use clean borosilicate glass bottles only.
- Prepare solvent volume to be consumed within 1 – 2 days.
- Use only HPLC-grade solvents filtered through 0.2  $\mu\text{m}$  filters.
- Label bottles correctly.
- Reduce risk of exposure to direct sun light by using brown bottles.

Preparing the Pump:

- Use fresh or different mobile phase (as required).
- Purge each channel with 5 mL/min for 10 min.
- Equilibrate your system (column and detector included) using the composition of your application for at least 15 min, until pressure and detector baseline signal are stable.

Daily Tasks:

- Replace water based mobile phases daily.
- Replace organic mobile phases at least every 2nd day.
- Check presence of seal wash solvent.
- Equilibrate the system with composition and flow rate of subsequent method.

Weekly Tasks:

- Change seal wash solvent (10 % isopropanol in water).
- Flush all channels with water at 2.5 – 3 mL/min for 10 min to remove salt deposits if buffer applications were used.
- Inspect solvent filters for dirt or blockages.

## Flushing Out Salt-Containing Solvents

- Replace the column with a union, replace the salt-containing solvent bottle with a new bottle of HPLC-grade water.
- Clean the bottle head assembly using lint-free wipes to minimize solvent carryover.
- Purge each channel for at least 10 min at 5 mL/min.
- Flush the entire system flow path with water for at least 10 min at 2 mL/min. During this step, switch the injection and the column selection valves between all positions at least 5 times.
- To minimize salt carryover, replace water with fresh solvent bottles.

## System Shutdown:

- Flush the column with the appropriate solvents and store it according to column manual instructions.
- Install a restriction capillary or union, and flush the system extensively with water, especially after using buffers.
- Flush and store the system in 50 % methanol or 50 % isopropanol in water, without additives.
- Power off the pump.

## Seal Wash (Usage Mandatory)

### 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:

The seal wash solvent is guided out to the waste on the right side of the module.

- PERIODIC operation, 0.5 min every 7 min when pump is switched on.
- Use 10 % isopropanol in water as seal wash solvent.
- Position wash solvent bottle above instrument.
- Exchange solvent weekly. Do not recycle seal wash solvent.

[www.agilent.com](http://www.agilent.com)

© Agilent Technologies Inc. 2021

Printed in Germany

Edition: 06/2021



Part No: G7132-90110 Rev. Rev. B.00

Document No: D0002930 Rev. B.00

