

**Ouick Reference Guide** 

# Maintenance Kits Make It Easy to Maintain Your LC Workflow Efficiency

Preventive maintenance helps stop problems like noisy baselines, short column life, and variable retention times. In fact, properly maintaining your LC system can lower your repair costs by 31%—and cut downtime by 2.4\* days per year.

Keep LC pumps, autosamplers, and detectors operating at their best with Agilent preventive maintenance kits. These contain the seals, frits, stators, fittings, connections, and other components you need to maximize LC instrument uptime.

Preventive maintenance kits are developed per module, so choosing a pump kit, autosampler kit, and detector lamp will cover most standard systems. We recommend scheduling pump and autosampler maintenance at a time when you are not using your instrument for analysis. For detectors, exchanging the lamp is sufficient.

\*Based on a five-year study of more than 38,000 service calls on Agilent and non-Agilent instruments, conducted by an independent actuarial firm. Read the complete white paper at <a href="https://www.agilent.com/chem/pm-whitepaper">www.agilent.com/chem/pm-whitepaper</a>.



## Need support?

Consult your system's user documentation for help with performing the required steps.

Find guidance at Agilent University: www.agilent.com/crosslab/university

Visit our resource page: www.agilent.com/chem/agilentresources

#### Pressed for time?

Several Agilent CrossLab service plans also cover preventive maintenance. Learn more: www.Agilent.com/en/support



# **Easy ordering information**

## Step 1

Identify which pump and autosampler your instrument uses. The module number appears at the lower-right corner on the front of each module.

#### Step 2

Use these tables to identify which kit corresponds to your pump and autosampler module.

## Step 3

Place your order at Agilent.com, or through your local Agilent representative.

### Preventive Maintenance Kits for Pumps

		G1310- 68741	G1310- 68742	G1312- 68741	G5611- 68741	G7104- 68741	G7120- 68741	G7161- 68730	G7161- 68740	G4204- 68741	G4204- 68742	G4220- 68741	G4220- 68742	G1310- 68730	G1310- 68731	G1312- 68730	G7120- 68708
System	Module	PM Kit	Add-on for Seal Wash	PM Kit	Add-on for Seal Wash	PM Kit	Add-on for Seal Wash	PM Kit	Add-on for Seal Wash	PM Kit	Universal Tool Kit						
1260 Infinity II	G7110B	•	•														•
	G7111A	•	•														•
	G7111B	•	•														•
	G7112B		•	•													•
	G7104C					•											•
	G5654A				•												•
1290 Infinity II	G7104A					•											•
	G7120A						•										•
	G7161A							•									•
	G7161B							•***	•								•
1290 Infinity	G4204A					• **				•	•						•
	G4220A						•**					•	•				•
	G4220B						•**					•	•				•
1260 Infinity	G1310B	•	•														•
	G1311B	•	•														•
	G1311C	•	•														•
	G1312B		•*	•													•
	G1312C		•*	•													•
	G5611A				•												•
1100/1200	G1310A													•	•		•
	G1311A													•	•		•
	G1312A															•	•

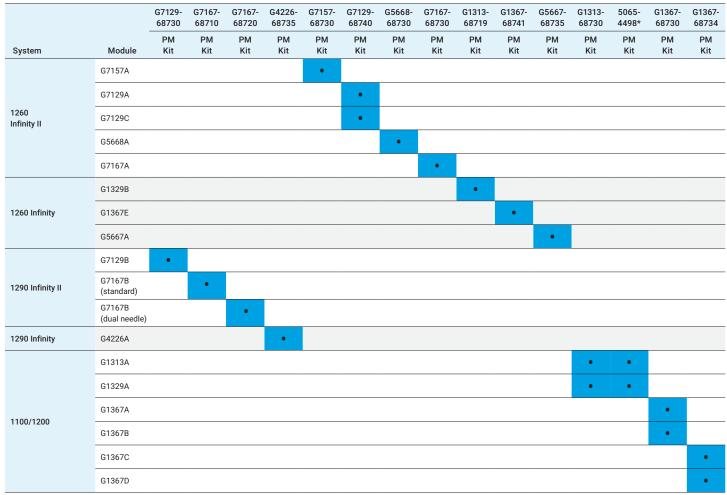
<sup>\*</sup>For G1312B and G1312C, two G1310-68742 add-ons are required.



<sup>\*\*</sup>Use this kit for pumps with upgraded "Easy Maintenance Pumpheads" (labeled as EM).

<sup>\*\*\*</sup>Use this kit when 50 mL/min pump heads are installed.

#### Preventive Maintenance Kits for Autosamplers



<sup>\*</sup>Extended preventive maintenance kit.

#### Preventive Maintenance Kits for 1120/1220 LC System

	G4280-68710	G4280-68730	G4280-68750	G4280-68770		
System	PM kit	PM kit	PM kit	PM kit		
1120 LC system with manual injector	•					
1120 LC system with automated injector		•				
1220 LC system with manual injector			•			
1220 LC system with automated injector				•		

# www.agilent.com/chem/store

DE44266.5072337963

This information is subject to change without notice.

