

Agilent 1290 Infinity

Binary Pump

Installation of the High Performance Jet Weaver

This technical note shows you how to install Jet Weaver for the 1290 Infinity Binary Pump.



Agilent Technologies

Introduction

The 380 μL Jet Weaver™ high performance mixer is optionally available for demanding applications, which use solvents in different channels (A vs. B), that differ strongly in their UV/Vis absorption, e.g. by using trifluoroacetic acid (TFA) as a modifier, which has a high absorbance. Solvent packages created by the pump may persist until the solvent reaches the detector flow cell. Absorption fluctuations can then show up as baseline noise, also referred to as mixing noise.

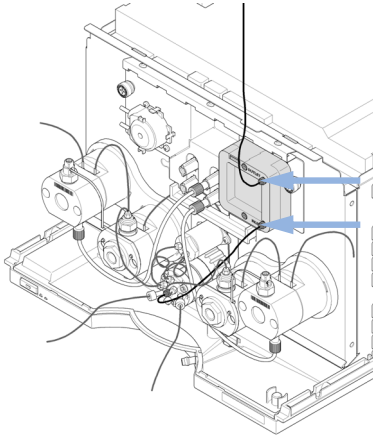
Applications like impurity quantitation or lowest level compound detection require minimizing this noise. The 380 μL Jet Weaver™ strongly improves mixing and therefore reduces baseline noise and improves sensitivity in detection. Patented Agilent microfluidic technology offers high mixing performance at a low internal volume of 380 μL (default mixer G4220-60006 has 35 μL or 100 μL).

Installation

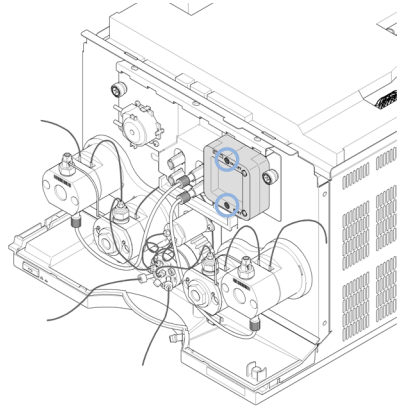
Tools required	p/n	Description
	8710-0510	¼ inch wrench
		3 mm hex key

Parts required	p/n	Description
	G4220-60012	Jet Weaver 380 µL
	G4220-87000	Capillary SST Valve to Jet Weaver 300 mm x 0.17 mm I.D.

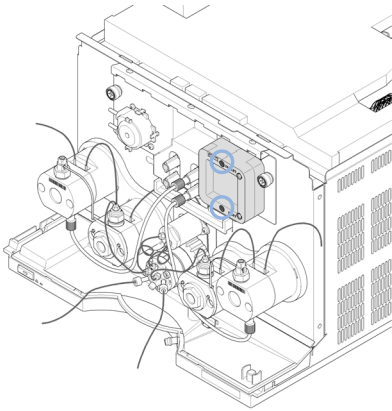
1 Remove capillary connections from the Jet Weaver.



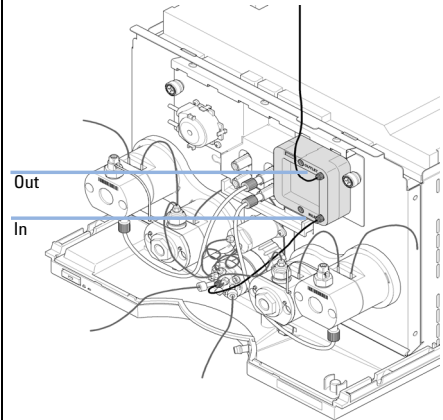
2 Remove the hex screws that fix the Jet Weaver to the pump housing.



3 Install new Jet Weaver by fixing it with the hex screws.



4 Reinstall the capillary connections.



The inlet at the bottom of the Jet Weaver is connected to the central port of the pump valve by a capillary (length 300 mm, 0.17 mm i.d.). The outlet at the top is connected to the autosampler.



G4220-90123

Part Number:
G4220-90123

Edition: 12/2011
Printed in Germany

© Agilent Technologies, Inc 2011

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