

## 11. Run test mix

Inject a test mixture to further measure system performance. The column test mixture used by Agilent to determine column quality is recommended. The Performance Summary Chromatogram included with each column can be easily duplicated if the same conditions and test mixture are used. Failure to duplicate the chromatogram for a new column indicates an installation, operation, or instrument problem. The problem must be corrected before proceeding with sample analysis.



## 12. Column use considerations

For maximum operating life, keep the column temperature below 100 °C when a column is installed but is not in use for analysis.

### Column storage

Seal the column ends with GC septa and return to the original box. Upon reinstallation, cut column ends to insure that no small pieces of septum have been left in the column.

### Chemical compatibility

Bonded and cross-linked stationary phases are not damaged by water or organic solvent injection. Inorganic acids (HCl, H<sub>2</sub>SO<sub>4</sub>, H<sub>3</sub>PO<sub>4</sub>, HNO<sub>3</sub>, etc.) and bases (KOH, NaOH, etc.) should not be injected into capillary columns. Rapid damage to the stationary phase will occur. If chemical damage does occur, removing the front ½ to 1 meter of the column will often restore column performance.

### Rinsing columns

*Do not solvent rinse the following non-bonded columns:* DX-1, DX-4, SE-30, SE-54, Carbowax 20M, Cyclodex-B, CycloSil-B, HP-20M, HP-88 and CAM, HP-101, HP-17, HP-Chiral β, CP-Sil 88, Select Silanes. CP Volamine, CP-TCEP, CP-Cyclodextrin, and CP-Chirasil Val.

All other Agilent J&W standard and cross-linked WCOT columns are solvent rinseable.

### Retention gaps

The stationary phase of non-bonded columns is easily disrupted during the injection process. Attach a 3 to 5 meter retention gap to the front of the column. This minimizes the amount of stationary phase damage, especially with on-column and splitless inlets.

## Temperature limits

Columns have both lower and upper temperature limits. Lower limits usually are at a phase change. Operation below this limit gives poor separation and peak shape problems, but will not damage the column.

Two upper limits are often given. The lower is the isothermal limit; the column can be held at this temperature for a prolonged time without harm. The higher one is a programming limit. The column can be heated to this limit for a short time (< 10 minutes).

Heating the column above the upper limits will significantly reduce column life. Set the GC oven maximum temperature at or below the column limit.

## The story behind Agilent J&W advanced GC columns

In 2000, Agilent, the inventor of fused silica GC tubing, acquired J&W Scientific, the creator of the first GC stationary phase made from cross-linked siloxane polymers. This partnership united the renowned DB and HP column families, creating Agilent J&W GC Columns. The innovative products that would follow include the Agilent J&W Ultra Inert, High Efficiency, LTM, and PAH GC columns. In 2010, Agilent acquired Varian, adding PLOT, Select, VF and CP-Sil to the most extensive and innovative GC column offering in the industry.

## The best low-bleed columns for sensitivity and performance

Column bleed can decrease spectral integrity, reduce uptime, and shorten column life. But Agilent J&W columns have the widest range of low-bleed and standard stationary phases featuring superior inertness and high upper temperature limits—especially for MS users performing trace level analyses.

## Better precision for better results

Agilent J&W columns adhere to tight retention factor (k) specifications, promoting consistent retention and separation. They also feature narrow retention indexes and a high number of theoretical plates per meter, ensuring narrow peaks and improving the resolution of closely eluting peaks.

## The industry's tightest quality control specifications

Agilent's stringent testing ensures reliable qualitative and quantitative results—even for your most challenging compounds. For example, we measure peak height ratios for both acids and bases to ensure top performance for the widest range of compounds. We also monitor peak symmetry and tailing for a broad scope of chemically active compounds.

As the world's leading provider of GC capillary columns, Agilent is uniquely positioned to offer you superior quality and unmatched service and support.

For additional column recommendations, chromatograms, and method parameters, go to [www.agilent.com/chem/myGCColumns](http://www.agilent.com/chem/myGCColumns)

## They may be small... but Agilent inlet supplies can have a big impact on your results

Agilent inlet supplies are engineered with the same reliability you expect from Agilent instruments. And they are designed to work with your new Agilent GC or GC/MS system. What's more, Agilent supplies are designed for fast, easy replacement, so you can get your system running again quickly.

Be sure to stock up on these essentials now—so they will always be there when you need them:

- Premium non-stick inlet septa and non-stick liner O-ring (5188-5365)
- Certified vials, caps and septa
- Ultra Inert injection port liners
- Pre-conditioned ferrules in ultra-clean packaging
- Metal injection-molded inlet Gold Seals (5188-5367)
- Gold standard autosampler syringes

**Learn more**  
[www.agilent.com/chem](http://www.agilent.com/chem)

**Buy online**  
[www.agilent.com/chem/store](http://www.agilent.com/chem/store)

**Find an Agilent Representative or Agilent Authorized Distributor**  
[www.agilent.com/chem/contactus](http://www.agilent.com/chem/contactus)

**U.S. and Canada**  
1-800-227-9770, [agilent\\_inquiries@agilent.com](mailto:agilent_inquiries@agilent.com)

**Europe**  
[info\\_agilent@agilent.com](mailto:info_agilent@agilent.com)

**Asia Pacific**  
[inquiry\\_lsca@agilent.com](mailto:inquiry_lsca@agilent.com)

**India**  
[india\\_lsca\\_marketing@agilent.com](mailto:india_lsca_marketing@agilent.com)

Agilent shall not be liable for errors contained herein or for incidental or consequential damages in connection with the furnishing, performance, or use of this material.

Information, descriptions, and specifications in this publication are subject to change without notice.

© Agilent Technologies, Inc., 2013  
Published in the USA, January 2, 2013  
Publication Number 830-0120



# AGILENT J&W CAPILLARY GC COLUMN INSTALLATION GUIDE

The Measure of Confidence

These instructions are suitable for the majority of the capillary columns that Agilent J&W manufactures. There may be more specific conditioning, care, and maintenance procedures for your new GC column (e.g., PLOT columns).

Be certain to read all the information that comes with your new GC column to ensure that the column performs to expectations.

A fused silica capillary GC column (**Figure 1**) consists of:

- An amber-brown polyimide exterior coating that protects the tubing from breakage.
- The fused silica tubing.
- A stationary phase that is evenly coated onto the inner wall of the tubing. Common phases are silicon-based polymers (polysiloxanes), polyethylene glycols, and solid adsorbents.

Maximize capillary GC column performance and lifetime by following these recommended guidelines for proper installation.

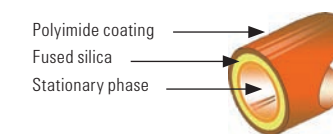
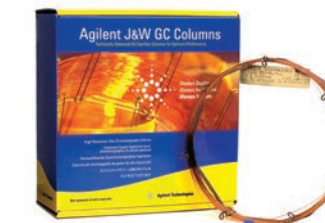


Figure 1. Capillary column



## Tools for capillary column installation

Column cutting tool such as a diamond-, carbide-, or sapphire-tipped pencil, or a ceramic cleaving wedge

Supply of an appropriate nonretained compound

Column test mixture (optional)

Electronic flowmeter (optional)

Electronic leak detector (optional)

## 1. Check traps, carrier gas, septum, and liner

Check gas purifiers for expiration and replace if necessary. Install a new septum in the inlet. If needed, clean or replace the inlet liner and gold-plated inlet seal, especially after injecting dirty samples or when analyzing active compounds. The liquid phase in the column is easily damaged by oxygen at temperatures above ambient. Use purifiers on the carrier gas lines to extend column lifetime and minimize background noise. An indicating filter that removes oxygen, moisture, and hydrocarbons such as the Agilent Gas Clean GC/MS filter is highly recommended for the carrier gas.



The Agilent Gas Clean filters can also be added to detectors gas lines to improve sensitivity and lifetime.

## 2. Place nut and ferrule on the column, carefully cut column end

**Warning:** Be careful! The oven and/or inlet may be hot enough to cause burns. If the inlet is hot, wear heat-resistant gloves to protect your hands.

**Warning:** Wear safety glasses to protect your eyes from flying particles while handling, cutting, or installing glass or fused silica capillary columns. Use care in handling these columns to prevent puncture wounds.

Place the column nut and ferrule over one end of the column. There is no front or back of the column; however, the posts of the column cage usually point towards the oven door.

Cut the end of the column after nut and ferrule placement. Hold the section of column to be cut against a finger. In one motion, scribe the outside of the column using a suitable cutting tool. Do not cut completely through the tubing. Grasp the column on each side of the scribe mark and bend away from the mark. Inspect the column end with a magnifier. Ensure the cut is at a right angle to the tubing wall and free of chips, burrs, or uneven areas (**Figure 2**). Recut if necessary.

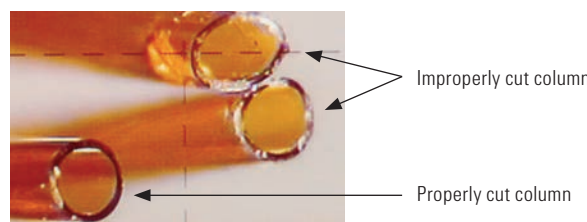


Figure 2. Cutting a fused silica capillary column.

## 3. Install column into the inlet

Place the column on the GC oven column hanger. Make sure the column tubing does not touch the sides of the oven.

Unwind enough column to obtain a smoothly curved section of tubing connected to the inlet. Avoid tight bends as this stresses the tubing and could cause breakage. Make sure that column tags, sharp edges, or other items do not rub against the column.

The optimal insertion distance of the column into the inlet depends on the inlet type. Consult the GC instruction manual for the proper insertion depth and technique.

With the column at its proper position, finger-tighten the column nut. Use a wrench to tighten an additional ½ turn. If the column can be moved in the fitting, tighten another ¼ turn. Failure to achieve a leak-free seal will cause rapid and permanent column damage. Do not move the column while tightening the nut.

Ferrules, especially those made of graphite/Vespel, will change shape slightly upon heating. If the column was installed while the inlet and detector were cool, retighten the fitting. It is also good practice to make sure the column nuts are tight after conditioning the column.

## 4. Turn on carrier gas

Adjust the head pressure to obtain a reasonable flow rate of carrier gas (**Table 1**). These values are recommended as starting points only. The actual head pressure will depend on the carrier gas velocity or flows set in Step 9.

Column length (m)	id (mm)						
	0.10	0.18	0.20	0.25	0.32	0.45	0.53
10	35-45	5-10					
12			10-15				
15				8-12	5-10		1-2
20	75-100	10-20					
25			20-30				
30				15-25	10-20	3-5	2-4
40		35-50					
50			30-60	15-25			
60				30-45	20-30	6-10	4-8
75						8-14	5-10
105			60-80				10-15

Table 1. Approximate column head pressure settings for He carrier gas (psig)

Place the free end of the column in a small vial of hexane. A steady stream of bubbles should be visible. If bubbles are not seen, check the carrier gas supply, flow controllers, etc. for proper settings and for leaks. Wipe off any residual solvent before continuing.

## Carrier gas selection

High purity helium and hydrogen are the preferred carrier gases for capillary columns; nitrogen is not recommended. Hydrogen is recommended for use with microbore GC columns (0.10 mm id and smaller). A gas purity of 99.995% or better is recommended with oxygen being the most important impurity to avoid (less than 1 ppm oxygen).

It is highly recommended to use gas purifiers such as the Agilent Gas Clean filter system, to ensure the highest quality gas while keeping gas lines clean and leak-free. Clean gases reduce risk of column damage, sensitivity loss, and instrument downtime.

**Warning:** Hydrogen forms explosive mixtures with air at concentrations of 4 to 10% hydrogen. Its high diffusivity minimizes this possibility, but the danger should not be discounted.

## 5. Install the column into the detector

Follow the installation precautions in Step 2 and 3 for the detector side while installing the column into the detector. Confirm all detector gas flows with an accurate flow-measuring device.

## 6. Inspect for leaks

Inspect the GC system for leaks before heating the column for the first time. An electronic leak detector is the most reliable way to check the inlet and detector fittings. Do not use Snoop. If use of a liquid is desired, try a 50/50 mixture of isopropanol/water.

## 7. Confirm carrier gas flow and proper column installation

Electronic pressure control (EPC) allows direct entry of carrier gas linear velocity or flow rate. It is critical that correct column dimensions are entered into the PC software or via the GC keypad for accurate velocity or flow values to be set.

Always consult the column performance summary sheet that accompanies the column for accurate inner diameter information.

Confirm carrier gas flow as described in Step 4 or by injecting a nonretained compound. Recommended nonretained compounds are in **Table 2**.

Detector	Compound
FID	Methane, butane
TCD	Methane, butane, argon, air
ECD	Methylene chloride <sup>2</sup> , SF <sub>6</sub> , CF <sub>2</sub> CL <sub>2</sub>
NPD	Acetonitrile <sup>2,3</sup>
PID	Ethylene, acetylene
MS	Methane, butane, argon, air

<sup>1</sup> Most of these compounds are significantly retained on PLOT columns.

<sup>2</sup> Do not inject liquid. Use a very dilute headspace injection.

<sup>3</sup> Acetonitrile is retained by most columns at temperatures below 100 °C. Heat the column to 100 °C or higher to set linear velocity.

Table 2. Nonretained compounds<sup>1</sup>

## Procedure

With the column temperature at 35 to 40 °C, rapidly inject 1 to 2 µL of a nonretained compound with the split/splitless inlet in the split mode. If using Megabore direct or cool on-column modes, dilute the nonretained compound so that the sample will not saturate the detector.

A very sharp and symmetrical peak should be obtained (**Figure 3**); a small amount of peak tailing may be observed with splitless injection. If no peak appears, there may not be any carrier gas flow. Check the regulators, gas supply and flow controllers for the proper settings. Make sure that the detector, recorder, and syringe are functioning properly. If the nonretained compound is tailing, there may be a leak in the inlet, poor column installation, or an excessively low split ratio. Reinstall the column and check the inlet for leaks. A nontailing peak is required before continuing.

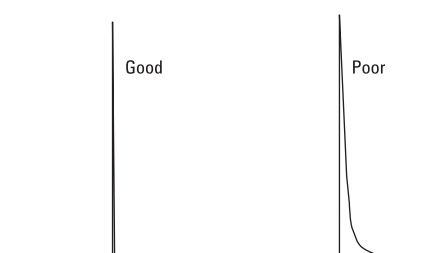


Figure 3. Methane peaks

## 8. Condition the column

**Caution:** Heating a column without carrier gas flow or while oxygen is present in the carrier gas stream will rapidly and permanently damage the column.

Purge the column with carrier gas for 15 minutes. Heat the column to its upper temperature limit or a temperature 10 to 20 °C above the highest operating temperature of the method, whichever is lower. Do not exceed the column upper limit or column damage will result.

After the column has reached the conditioning temperature, observe the baseline. It will rise for 5 to 30 minutes, then drop for another 30 to 90 minutes. A flat baseline should be obtained 1 to 3 hours after reaching the conditioning temperature. If the baseline does not stabilize after 2 to 3 hours or does not remain constant, stop the conditioning process.\* An unstable baseline can be caused by a leak in the carrier gas line or inlet area, or by system contamination. Fix either problem before continuing.

Polar stationary phases and thick films usually require longer times to stabilize than nonpolar phases and thinner films. PLOT columns require special conditioning procedures. Refer to the column information sheet for the appropriate procedure.

Some detectors, such as ECDs and MSDs, may stabilize faster if the column is not connected to the detector during column conditioning.

If the column is conditioned with the detector end disconnected, a small portion of the column end may be damaged. Remove 10 to 20 cm of the exposed column end before connecting the column to the detector.

\* Thick-film and PLOT columns may take longer.

## 9. Accurately set the carrier gas velocity

For capillary columns, the average linear velocity (µ) is a better and more meaningful measure of the carrier gas than the volumetric flow rate. The carrier gas linear velocity directly influences the retention time and efficiency.

## GCs without EPC

Carrier gas velocity changes with oven temperature as carrier gas viscosity changes. Always set linear velocity at the same temperature for a given analysis (often the initial oven temperature). Inject 1 to 2 µL of the appropriate nonretained compound and calculate the linear velocity using the retention time of the peak and the following

equation. Adjust the column head pressure until the desired average linear velocity is obtained.

$$\mu = L/t_r$$

Where  
 $\mu$  = Average linear velocity (cm/sec)  
 $L$  = Column length (cm)  
 $t_r$  = Retention time nonretained peak (sec)

## Recommended average linear velocities

Helium:	30 to 40 cm/sec
Hydrogen:	50 to 80 cm/sec

## GCs with EPC

The linear velocity can be entered into the PC software or the GC keypad and remains constant. The correct column dimensions must be entered into the PC software or the GC keypad for the system to accurately set the linear velocity. Refer to the inner diameter of the column which is included on the column performance summary sheet (this is a nominal value).

## 10. Bleed test

After the column is conditioned, run a blank (no injection) chromatogram. Start at 40 to 50 °C, ramp at 10 to 20 °C/min and hold for 10 to 15 minutes at the conditioning temperature. Save this background trace for future comparisons. See **Figure 4**.

There should be no peaks in the blank chromatogram. Peaks indicate a contamination problem, usually in the inlet area. As a column degrades with normal usage, the magnitude of the baseline rise will increase. If the baseline rise occurs at a much lower temperature than previously obtained, the column and/or GC is most likely contaminated or damaged.

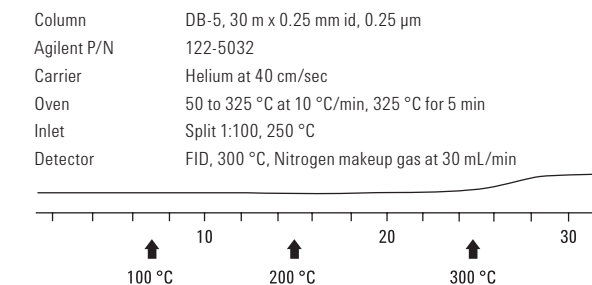


Figure 4. Bleed profile