



## GC Column Installation Quick Reference Guide - Inlets

Inlet	Diagram	Procedure
Split/Splitless		<p>Place a septum over the column, then the column nut and ferrule. Trim the end of the column with a column cutter.</p> <p>Pull the column back so that 4-6 mm of column is extending past the end of the ferrule.</p> <p>Thread the column nut and column into the inlet and tighten slightly past where the column grabs – retighten after heating.</p>
Purged Packed		<p>Place a septum over the column, then the column nut and ferrule. Trim the end of the column with a column cutter.</p> <p>Pull the column back so that 1-2 mm of column is extending past the end of the ferrule.</p> <p>Thread the column nut and column into the inlet and tighten slightly past where the column grabs – retighten after heating.</p>
Multimode		<p><b>NOTE:</b> Make sure the column adapter nut on the inlet base is <b>fully threaded on</b> and <b>spinning freely – Collar Up!</b></p> <div style="text-align: center;"> </div> <p><b>Tighten with two wrenches</b> - 1/4" and 5/16" To prevent damage the inlet threads.</p>
Cool On Column		<p>Insert the column all the way into the inlet until you feel the spring tension – do not withdraw.</p> <p><b>The column cut is critical.</b></p> <p><b>Tighten with two wrenches</b> - 1/4" and 5/16" to avoid damaging the inlet.</p>
PTV		<p>There should be 17mm of column above the graphpak ferrule – the graphpak ferrule should be installed with the graphite end towards the inlet base. The column nut is slotted. Use a 5 mm wrench to tighten the fitting.</p>
Volatiles Interface		<p>There is a longer column nut for the VI so that you don't have to remove the inlet block.</p> <p>Part Number - G3504-20504</p>



## GC Column Installation Quick Reference Guide - Detectors

Detector	Diagram	Procedure
FID/NPD		<p>Place a septum over the column, then the column nut and ferrule. Trim the end of the column with a column cutter.</p> <p>Thread the column nut and column into the detector base. Insert the column all the way into the detector jet until it stops, then withdraw 1-2 mm before tightening the nut.</p> <p>The dimensions shown of 48 and 68 mm are provided for very narrow columns (&lt;100um) which may go all the way through the jet.</p>
TCD – Using the standard Low Leakage Ferrules		<p>Place a 1/8" Swagelok nut and the Back and Front ferrules over the column. (Be sure to select the front ferrule size to match the column) Trim the end of the column with a column cutter.</p> <p>Thread the column nut and column into the detector base. Insert the column all the way into the detector jet until it stops, then withdraw 1-2 mm before tightening the nut. Re-tighten after heating.</p>
TCD – Original Design Column Adapter		<p>Install the column adapter to the 1/8" Swagelok TCD base using a graphite/vespel ferrule.</p> <p>Thread the column nut and column into the adapter. Insert the column all the way into the detector jet until it stops, then withdraw 1-2 mm before tightening the nut.</p>
uECD		<p>Be sure that the Makeup gas adapter is installed with the 19 mm dimension shown.</p> <p>Thread the column nut with column into the make-up gas adapter. Insert the column all the way into the mixing liner until it stops, then withdraw 1-2 mm - Use 2 wrenches to tighten the nut. Re-tighten after heating.</p> <p>The 70 mm dimension shown is provided for very narrow columns (&lt;100um) which may go all the way through the mixing liner.</p>
FPD		<p>Install the column into the Column Measuring Tool Provided. Trim the column such that 1 mm extends from the end of the tool. Tighten the column nut and mark the position of the column at the back of the nut. Install into the detector base.</p>