

Finger-Tight Column Nut with Internal Threads Use Instructions

PerkinElmer's Finger-Tight Column Nut with Internal Threads (PN: N6400721) is a fitting designed for easy toolless column installation. The fitting includes a disk spring that loads as you tighten the fitting. Using Graphite-Vespel ferrules with the Finger-Tight nut is recommended. This pairing ensures a tight seal that requires no retightening, even after successive thermal cycles.

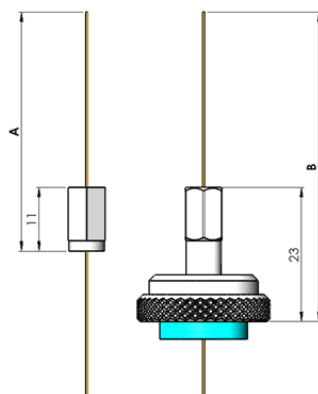
Compatible Injectors and Detectors

The Finger Tight Column Nut with Internal Threads is compatible with the inlets of the GC 2400 PerkinElmer capillary PSS injector (N6400051) and Mass Spectrometer MS 2400 (N6480023).

Recommended PerkinElmer Ferrules

Part Number	Description	ID	Material	Holes
04972392	FERRULE VG 1/16 STD 2 HOLE 0.4MMID 10PK	0.4	85% Graphite / 15% Vespel	2
09920104	FERRULE STD VG 1/16 0.4MM ID 10PK	0.4	85% Graphite / 15% Vespel	1
N9307245	FERRULE STD GRAPH 1/16 0.4MM ID 10PK	0.4	Graphite	1
L1003027	FERRULE STD VG 1/16 0.5MM ID EACH	0.5	85% Graphite / 15% Vespel	1
04972433	FERRULE STD GRAPH 1/16 0.5MM ID EACH	0.5	Graphite	1
09200788	FERRULE VG 1/16 STD 2HOLE 0.5MMID 10PK	0.5	85% Graphite / 15% Vespel	2
09903700	FERRULE STD GRAPH 1/16 0.5MM ID 10PK	0.5	Graphite	1
09920105	FERRULE STD VG 1/16 0.5MM ID 10PK	0.5	85% Graphite / 15% Vespel	1
09920107	FERRULE STD VG 1/16 0.8MM ID 10PK	0.8	85% Graphite / 15% Vespel	1
09920141	FERRULE STD GRAPH 1/16 0.8MM ID 10PK	0.8	Graphite	1
N6710240	FERRULE STD VG 1/16 0.9MM ID 10PK	0.9	85% Graphite / 15% Vespel	1
N9307244	FERRULE STD GRAPH 1/16 1.0MM ID 10PK	1.0	Graphite	1
04972263	FERRULE STD VG 1/16 1.6MM ID EACH	1.6	85% Graphite / 15% Vespel	1
09920127	FERRULE STD VG 1/16 1.6MM ID 10PK	1.6	85% Graphite / 15% Vespel	1
09920511	FERRULE VG NO-HOLE 1/16 STD 10PK	0.0	85% Graphite / 15% Vespel	0

Column Lengths of Finger-Tight Fittings



Column Length (mm) from Back of the Fitting

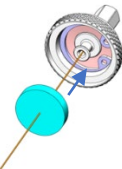
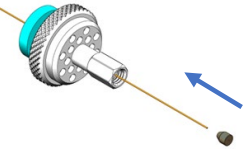
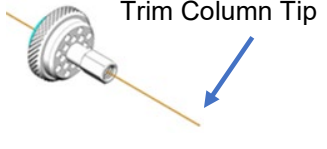
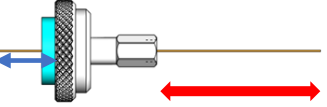
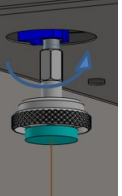
Device	09903392 [A]	N6400481 [B]
PSS Injector	38-44	50-56
MS 2400	See MS 2400 Instructions	See MS 2400 Instructions



PerkinElmer U.S. LLC
710 Bridgeport Avenue
Shelton, CT 06484-4794, U.S.A

Finger Tight Use Instructions

To install a column or detector with a Finger Tight Fitting

<ol style="list-style-type: none"> 1. Push a septum (N9306219 or similar) onto the GC column from the column tip. 2. Push the column tip into the Finger Tight fitting, so that the column fits snugly into the fitting. 	
<ol style="list-style-type: none"> 3. Install a long ferrule (minimum 3.7mm height 2-hole ferrules acceptable) onto the column, making sure that the tapered end of the ferrule points into the fitting. <p>Note: For more information, see Recommended Ferrule Types on the reverse side.</p>	
<ol style="list-style-type: none"> 4. After inserting the column through the ferrule, trim about 1 centimeter from the tip of the column for a clean, debris free column. Use the Wafer-Ceramic Cutter (N9301376) to scribe the column end. Then break off the tubing at the score mark so that the break is clean and square. 	 <p>Trim Column Tip</p>
<ol style="list-style-type: none"> 5. Adjust the location of the septum along the column (blue arrow) to place the tip of the column at the appropriate length (red arrow) for the injector or detector being used. <p>Note: For more information, see Column Lengths of Finger-Tight Fittings on reverse side.</p>	
<ol style="list-style-type: none"> 6. Insert the tip of the column into the injector or detector. 7. Holding the septa against the septa stop, tighten the fitting. <p>Note: When the fitting begins to crimp the ferrule, tighten the fitting another one quarter turn. Confirm that the ferrule is fully crimped onto the GC column by gently pulling on the GC column. It should not slip.</p>	

Removing a Column with a Finger-Tight Fitting

To remove a column installed with a Finger-Tight fitting:

1. Grip the Finger-Tight fitting and back it out of the GC 2400 PSS injector or MS2400 inlet.
2. Pull Graphite-Vespel ferrule off the tip of the fitting.

Note: Graphite ferrules may require a pointed tool to dislodge them from the fitting ferrule cup.