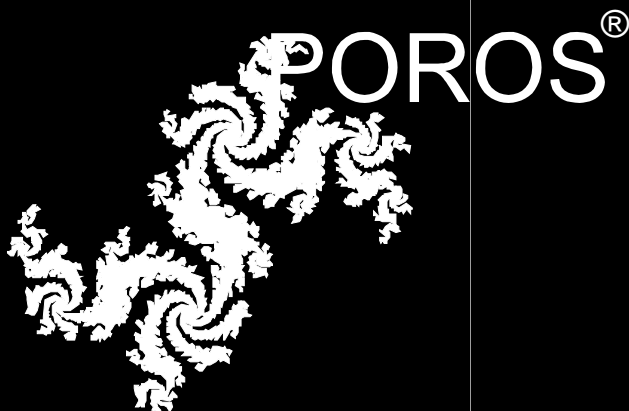


# Inlet Frit Replacements PEEK Column



**PE Biosystems  
Headquarters  
850 Lincoln Centre Drive  
Foster City, CA 94404  
Tel: (650) 638-5800  
Tel: (800) 345-5224  
[www.pebiosystems.com](http://www.pebiosystems.com)**

**PerSeptive Biosystems  
Center of Excellence  
500 Old Connecticut Path  
Framingham, MA 01701  
Tel: (508) 383-7700  
Tel: (800) 899-5858  
[www.pbio.com](http://www.pbio.com)**

For international office locations, please call the headquarters or refer to our web site at [www.pebiosystems.com](http://www.pebiosystems.com).



**PerSeptive Biosystems**

PerSeptive Biosystems, POROS, Perfusion Chromatography, and the fractal icon are registered trademarks of PerSeptive Biosystems, Inc.

Part Number 8-0034-40-0993, Rev. 3, 9/95 ©1995 Printed In USA



Printed on recycled paper.

To replace the inlet frit:

1. Seal the outlet end of the column with an end plug.
2. Carefully unscrew the inlet end fitting from the column and remove it by hand. If necessary, grip the column tube with a rubber glove while you remove the end fitting with a wrench.
3. Remove the frit from the end fitting (or the end of the column) and flush out excess packing material from the fitting.
4. Hold the end fitting with the opening facing up and drop the new frit into the fitting.
5. Keeping the end fitting in the same position, turn the column upside down and screw it into the fitting until it is finger-tight. Then gently tighten it with a wrench.
6. Backflush the column at low flow until all the air is removed.
7. Check for particles in the effluent to ensure that the frit is properly seated.
8. If the column leaks, tighten the end fitting slightly.
9. After buffer conditioning, check column pressure and efficiency.