



G-Biosciences ♦ 1-800-628-7730 ♦ 1-314-991-6034 ♦ technical@GBiosciences.com

A Geno Technology, Inc. (USA) brand name

G-Trap™ FPLC Columns

Datasheet for the 1 & 5ml FPLC Columns

(Cat. # 786-1290, 786-1291)



think proteins! think G-Biosciences www.GBiosciences.com

ITEM(S) SUPPLIED

Cat. #	Description	Size
786-1290	G-Trap™ 1ml FPLC Column	1
786-1291	G-Trap™ 5ml FPLC Column	1

SPECIFICATIONS

- **Column material:** polypropylene (PP)
- **Frit material:** Ultra high molecular weight polyethylene (UHMW-PE)
- **Maximum pressure:** 3 bar (44 psi)

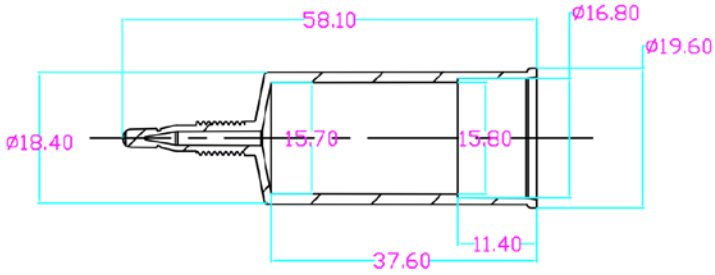
COMPONENTS AND DIMENSIONS

- 1 Column tube
- 1 Sealing plug
- 1 Sealing sleeve
- 1 Top stop plug & 1 bottom stop plug
- 1 Top (Large) frit
- 1 Bottom (Small) frit (inserted in column)

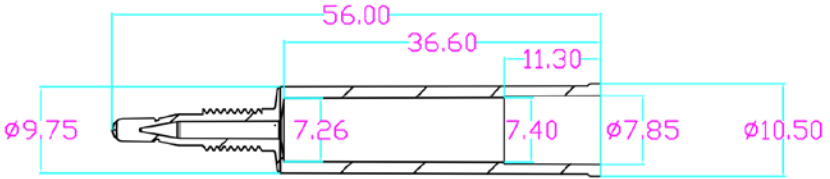


1. Place bottom frit into the column and slide to the bottom.
NOTE: The bottom frit may already be present in the column
2. Fill column with 1 or 5ml resin as per resin manufacturer's guidelines.
NOTE: Avoid air bubbles.
3. Apply top frit and position above the resin with no gap.
4. Add the sealing plug and push firmly into column.
5. Add the sealing sleeve and seal the column with the Top Stop Plug.

Column Tubes

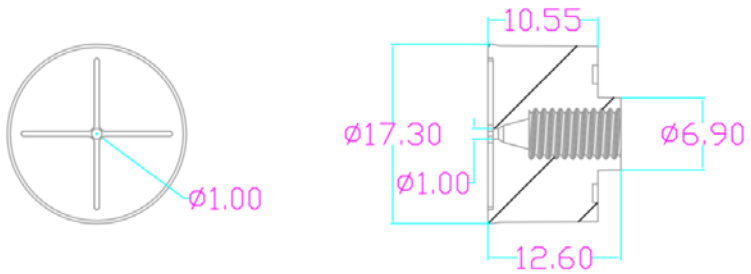


5-mL column tube

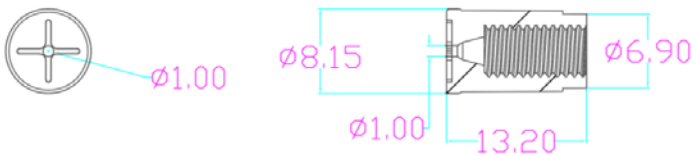


1-mL column tube

Sealing Plugs

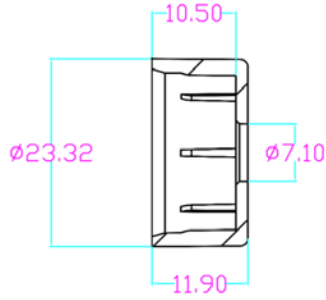


Sealing plug for 5-mL column, showing bottom view (left) and side view (right)

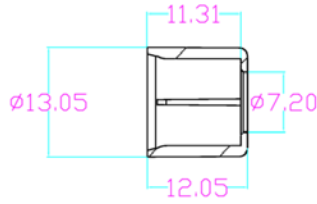


Sealing plug for 1-mL column, showing bottom view (left) and side view (right)

Sealing Sleeves

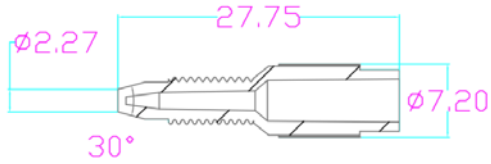


Sealing sleeve for 5-mL column

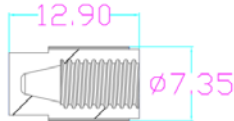


Sealing sleeve for 1-mL column

Top and Bottom Stop Plugs



Top stop plug (male finger-tight stop plug)



Bottom stop plug (female finger-tight stop plug)

RELATED PRODUCTS

For other related products, visit our website at www.GBiosciences.com or contact us.

Last saved: 8/23/2017 IA



www.GBiosciences.com