

EconoFit Empty Columns, 1 and 5 ml

Catalog #	Description
12009232	EconoFit Empty Columns , 5 x 1 ml
12009233	EconoFit Empty Columns , 5 x 5 ml

For research purposes only.

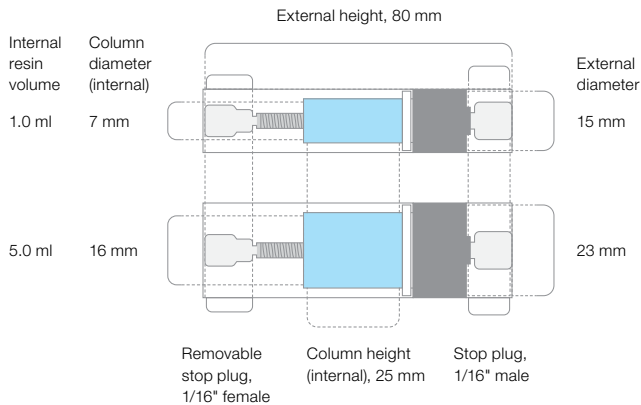
Product Information

EconoFit Columns are disposable, easy-to-use, prepacked chromatographic columns supplied ready for use in convenient 1 and 5 ml sizes. They are quickly connected to liquid chromatography systems using 10-32 fittings. Columns are available for a variety of chromatographic techniques, including desalting (size exclusion), ion exchange, affinity, mixed-mode, and hydrophobic interaction chromatography. See Table 1 for specifications.

Table 1. EconoFit Empty Column specifications.

Property	Description
Size	1 and 5 ml bed volumes
Dimensions	1 ml: 80 mm length x 15 mm external diameter 5 ml: 80 mm length x 23 mm external diameter
Recommended flow rate (depending on resin type)	1 ml: 1–2 ml/min (240–480 cm/hr) 5 ml: 5–10 ml/min (240–480 cm/hr)
Maximum flow rate	1 ml: 6 ml/min (1,440 cm/hr) 5 ml: 20 ml/min (963 cm/hr)
Fittings	10-32 (1/16"), female inlet and male outlet
Column material	Polypropylene
Frit material	High-density polyethylene
Storage conditions	20% ethanol
Autoclavability	Not autoclavable

BIO-RAD is a trademark of Bio-Rad Laboratories, Inc.
All trademarks used herein are the property of their respective owner.



Schematic of EconoFit Columns.

Chemical Compatibility

The EconoFit Column hardware is resistant to the majority of commonly used buffers and reagents used in the purification of biomolecules. These include, but are not limited to, the following:

- pH range 1–14 (short-term exposure at extreme pHs)
- 2% benzyl alcohol
- 20% ethanol
- Up to 1 M NaOH
- Up to 1 M nonoxidizing mineral acids
- Detergents
- Ethylene glycol
- 6 M guanidine HCl
- 8 M urea
- Nonhalogenated organic solvents (for example, acetone, acetonitrile)
- Reducing agents (such as DTT) and chelating agents (such as EDTA)

For full details of all prepacked EconoFit Columns, refer to the EconoFit Chromatography Columns Product Insert (10000102129).

Visit bio-rad.com/ResinsandColumns for a complete listing of items in the EconoFit Column product line.