

Agilent Packed GC Column user manual



Packaging

Agilent J&W Packed GC Columns are packaged in a special bag, which prevents moisture and oxygen from entering the column. This ensures the cleanliness of the packed GC column.

IMPORTANT: Please make sure to remove yellow plastic caps from column legs before installing the column into the GC inlet and detector.

Preconditioning

Heat the column 20 °C below the maximum temperature.

Exceptions are:

Molecular Sieve 5A and Molecular Sieve 13X Columns.

Heat the column from 30 °C to 300 °C at 0.5 °C/min.

Use a purified gas with a minimum flow of 20 mL/min.



Agilent Technologies

Gas purification

The use of Agilent Gas Clean filters with indicators is strongly recommended. Agilent Gas Clean filters are designed to prevent all types of contamination from causing damage to the column, loss of sensitivity and unnecessary instrument downtime. Agilent's unique, high sensitivity indicators warn you if contaminants are present and when it is time to replace the filters.

Optimum flow

Approximate flow as a relation of carrier gas and inside diameter

id	Nitrogen (mL/Min)	Helium, Hydrogen (mL/Min)
2 mm	15	20
3 mm	20	25
4 mm	25	30

Ordering information

Agilent Gas Clean Filters

Description	Part No.
Agilent Gas Clean CO ₂ Filter	CP17969
Agilent Gas Clean Oxygen Filter	CP17970
Agilent Gas Clean Moisture Filter	CP17971
Agilent Gas Clean Process Moisture Filter	CP17971P
Agilent Gas Clean Charcoal Filter	CP17972
Agilent Gas Clean GC/MS Filter	CP17973

www.agilent.com/chem

Agilent shall not be liable for errors contained herein or for incidental or consequential damages in connection with the furnishing, performance, or use of this material.

Information, descriptions, and specifications in this publication are subject to change without notice.

© Agilent Technologies, Inc. 2015
Published in The Netherlands,
February 11, 2015
Part Number 5991-5578EN

