

FIND

YOUR IDEAL MATCH

The Measure of Confidence



Agilent Parts Finder Tool and GC Column Selection Tool



Satisfaction GUARANTEED!

Quickly locate and order the Agilent supplies you need with our Parts Finder Tool

This intuitive wizard features an image-based search and inventory for quick identification of supplies for your Agilent GC instruments.

- Select GC supplies in seconds
- Quickly locate hard to find part numbers and build a list of favorites
- Print your parts list or save it to a file

Try it now at agilent.com/chem/partsfinder



Agilent Technologies



Which Agilent J&W GC column is right for you? Find out in seconds with our GC Column Selection Tool

Let the chromatography experts help you select the right GC column – right from your computer or Smartphone. And to make it even easier, we give you three ways to search:

1. By method
2. By column phase, length, id, and film thickness
3. By Agilent or non-Agilent part number

Get started now at agilent.com/chem/SelectGC

SPECIAL SAVINGS

With the purchase of a qualifying Agilent J&W column or Agilent supply. Limited time offer.

Shop now at agilent.com/chem/specialoffers

This information is subject to change without notice.

©Agilent Technologies, Inc., 2013

Printed in the USA, May 5, 2013

5991-2321EN

