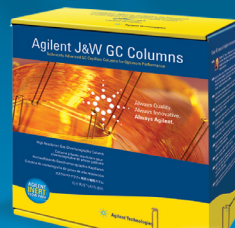


IMPROVED AGILENT J&W HP-INNOWAX GC COLUMNS

ENHANCE YOUR POLAR PEAK SHAPE ACCURATE. SENSITIVE. REPEATABLE.



Now your analysis of polar compounds can be even more worry free

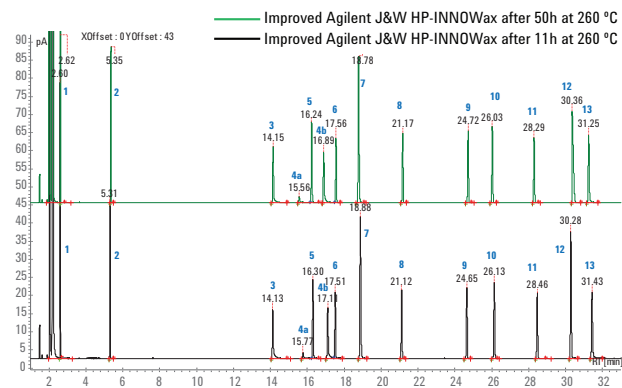
For increasingly active polar analytes, you cannot afford adsorption caused by flow path activity. The Agilent J&W Ultra Inert GC column family pushes industry standards for consistent column inertness and exceptionally low column bleed. The innovative processes, employed in the manufacture of Agilent J&W DB-Wax Ultra Inert columns, are now being applied to the production process for the HP-INNOWax GC Columns.

As a result, the Improved Agilent J&W HP-INNOWax GC columns now deliver:

- Excellent peak shape performance for active polar compounds
- Better inertness longevity performance
- Improved column-to-column inertness reproducibility and retention time stability
- Extended inertness lifetime that withstands repeated cycling to the upper temperature limits of the column

Other key performance parameters—such as selectivity, theoretical plates, and retention indices—remain unchanged. This makes the transition to the new improved columns seamless.

Excellent peak shape for Grob test mixture



HP-INNOWax GC columns show outstanding performance for this modified Grob mixture even after exposure to the upper temperature limit for 50 hours.

Peak no.	Compound	Peak no.	Compound
1	Decane	8	nC11-FAME
2	Dodecane	9	nC12-FAME
3	Decanal	10	2,6-Dimethylaniline
4	2,3-Butanediol**	11	2,6-Dimethylphenol
5	1-Octanol	12	2-Ethylhexanoic acid
6	C10-FAME	13	Ethyl maltol
7	Dicyclohexylamine		


** 2,3-Butanediol is present at 2 isomers: RR/SS and meso isomers, respectively.



Agilent Technologies

IMPROVED AGILENT J&W HP-INNOWAX GC COLUMNS ARE PART OF THE AGILENT ULTRA INERT GC FLOW PATH

By minimizing activity along every step of the GC and GC/MS flow path, Agilent Inert Flow Path solutions improve system performance, ensure better results, and allow you to process more samples without unplanned maintenance and recalibration. So you won't miss a thing in your GC analysis.



Ultra Inert liners



Inert Flow Path Split/Splitless inlet



Ultra Inert gold seals



Inert MS source



IDP-3 pump

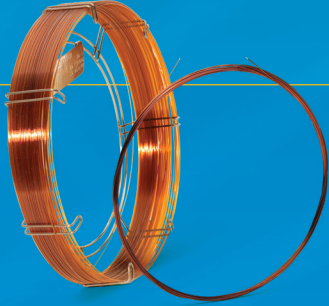





UltiMetal Plus Flexible Metal ferrules



Inert Capillary Flow Technology devices, including UltiMetal Plus 3-way splitter



Improved Agilent J&W HP-INNOWax GC column and Ultimate Plus deactivated fused silica tubing



Gas Clean purifier

Ordering Guide

ID (mm)	Length(m)	Film(μm)	Temp Limits (°C)	7 in Cage	5 in Cage	7890/6890 LTM II module
0.18	20	0.18	40 to 260/270	19091N-577i	19091N-577iE	
0.20	25	0.20	40 to 260/270	19091N-102i		
	50	0.20	40 to 260/270	19091N-105i		
		0.40	40 to 260/270	19091N-205i		
0.25	15	0.25	40 to 260/270	19091N-131i		
		0.50	40 to 260/270	19091N-231i		
	30	0.15	40 to 260/270	19091N-033i		
		0.25	40 to 260/270	19091N-133i	19091N-133iE	19091N-133iLTM
		0.50	40 to 260/270	19091N-233i	19091N-233iE	
	60	0.25	40 to 260/270	19091N-136i	19091N-136iE	
		0.50	40 to 260/270	19091N-236i		
0.32	15	0.25	40 to 260/270	19091N-111i		
	30	0.15	40 to 260/270	19091N-013i		
		0.25	40 to 260/270	19091N-113i	19091N-113iE	
		0.50	40 to 260/270	19091N-213i	19091N-213iE	
	60	0.25	40 to 260/270	19091N-116i		
		0.50	40 to 260/270	19091N-216i	19091N-216iE	
0.53	15	1.00	40 to 240/250	19095N-121i		
	30	1.00	40 to 240/250	19095N-123i	19095N-123iE	
	60	1.00	40 to 240/250	19095N-126i		

Learn more about analyzing polar compounds with utmost confidence
www.agilent.com/chem/hp-innowax

For Research Use Only.
Not for use in diagnostic procedures.
This information is subject to change without notice.

© Agilent Technologies, Inc., 2016
Printed in the USA, December 6, 2016
5991-7635EN

