

Selecting the Right WAX Column

Agilent J&W WAX GC columns deliver excellent inertness with greater sensitivity than competitive WAX columns. With the increasing number of available Agilent J&W WAX columns, it can be challenging deciding which one is best suited for your needs.

This poster will help you select the right WAX column for optimal performance, based on factors such as application analyte mix and the type of detector being used.

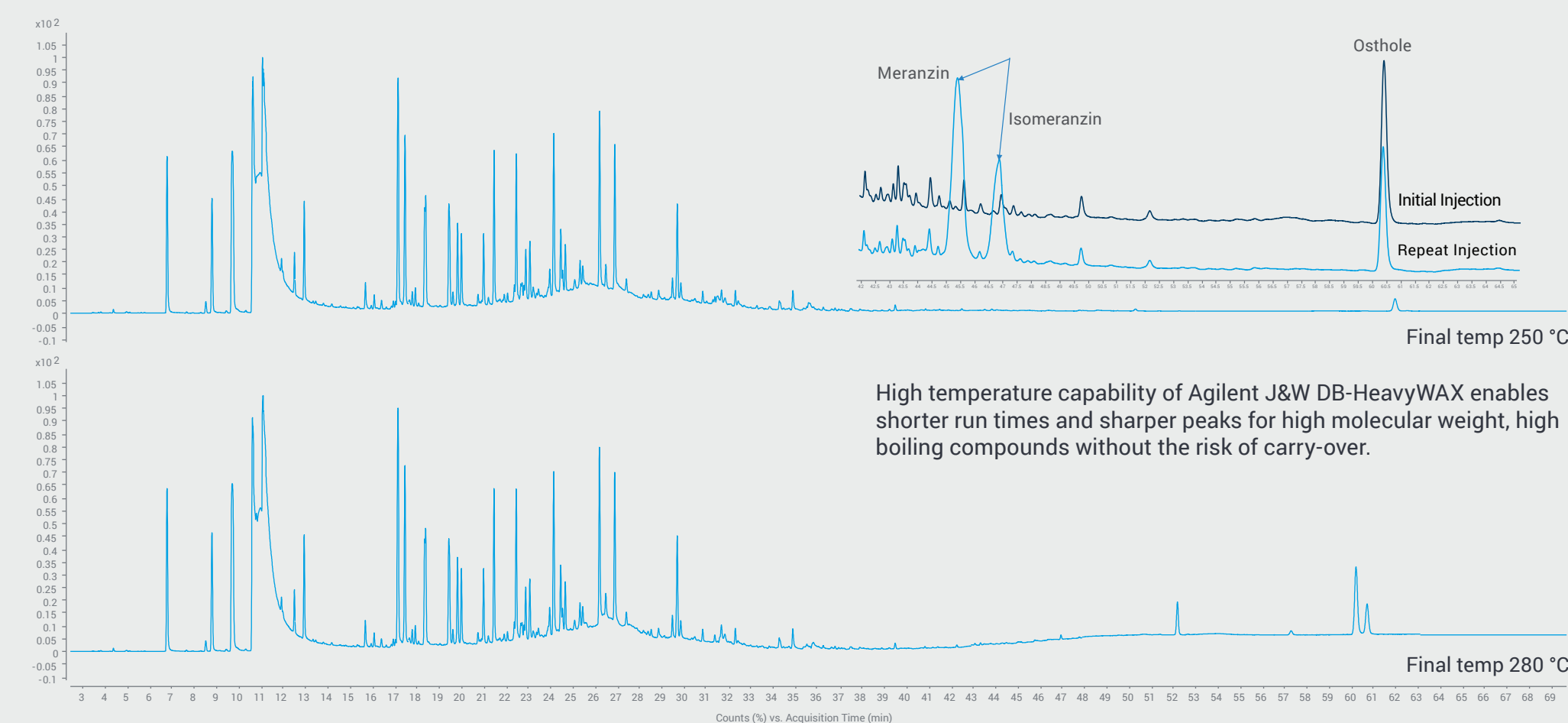


DB-HeavyWAX for high molecular weight analytes

- Higher max temp limit: 280 °C isothermal and 290 °C programmed
- Shorter run times

www.agilent.com/chem/db-heavywax

Analysis of Cold Pressed Pink Grapefruit Essential Oil



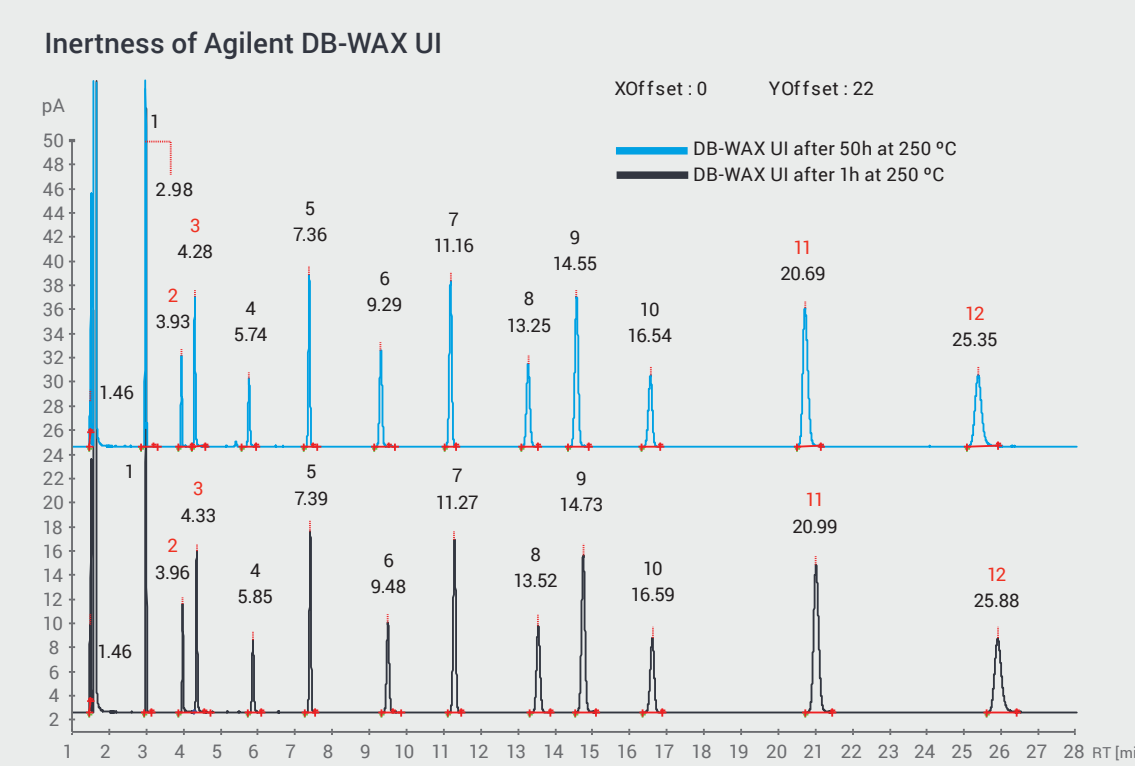
DB-WAX Ultra Inert for polar compounds



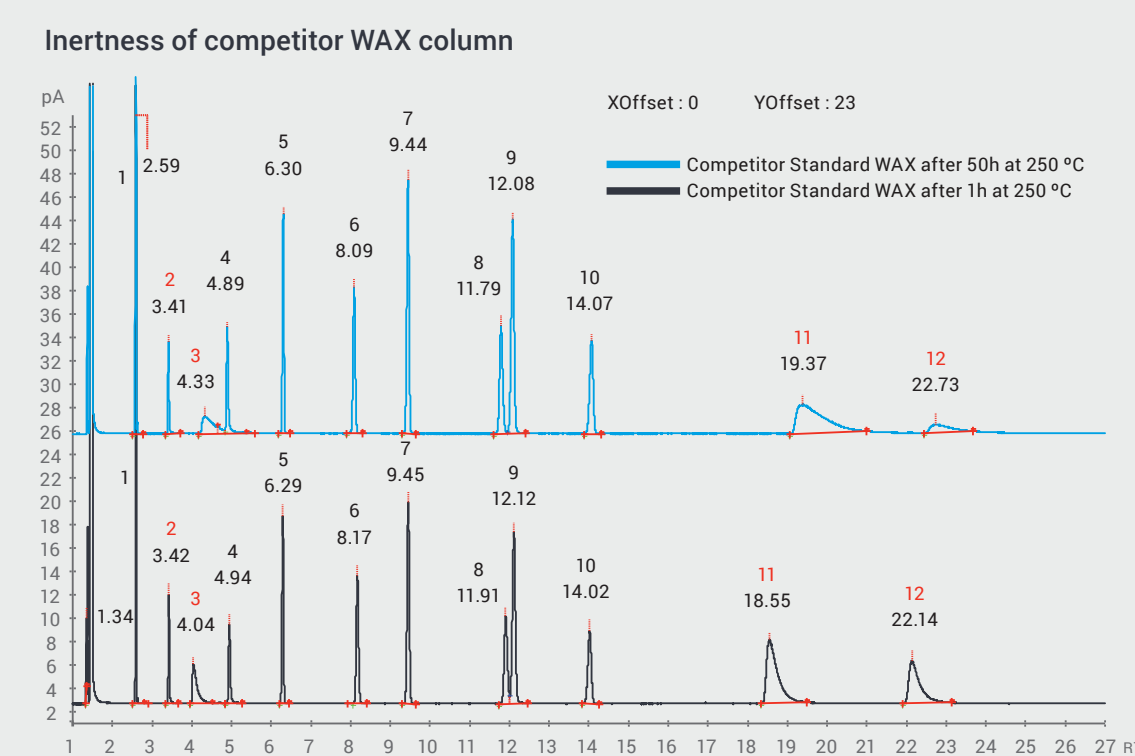
- Withstands repeated temperature cycling to upper limits without sacrificing peak shape performance
 - Identical selectivity to DB-WAX
- www.agilent.com/chem/dbwaxui

1. 2-Nonanone
2. Decanal
3. Propionic acid
4. Ethylene glycol
5. Heptadecane
6. Aniline
7. Methyl dodecanoate
8. 2-Chlorophenol
9. 1-Undecanol
10. Nonadecane
11. 2-Ethylhexanoic acid
12. Ethyl maltol

Injector: 250 °C, Split 1:75
Injection volume: 1 µL
Flow rate: 1.1 mL/min, H2 gas
Oven program: 130 °C isothermal
Detector: FID @ 260 °C



Even after 50 hours of exposure at the upper temperature range, the DB-WAX UI GC column demonstrated superior inertness and stability with minimal peak shifting.



DB-FATWAX Ultra Inert for FAMES and free fatty acids

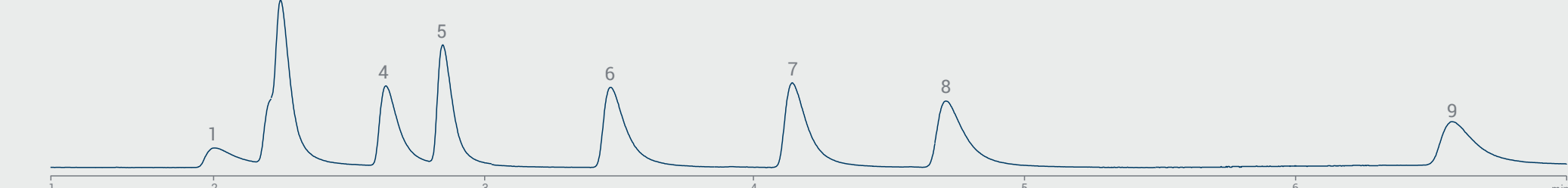


- Single column analysis
 - Solvent rinseable and tolerates aqueous injections
- www.agilent.com/chem/fatwax-ui

Superior inertness, unique selectivity, and thermal robustness for separation of FAMES and underivatized free fatty acids.

Competitor WAX column

30m x 0.25 mm, 0.25 µm



Agilent DB-FATWAX Ultra Inert

30m x 0.25 mm, 0.25 µm

