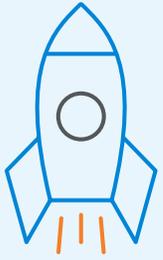


# Built for Ultrashort-Chain PFAS. Compatible for C1 to C18.

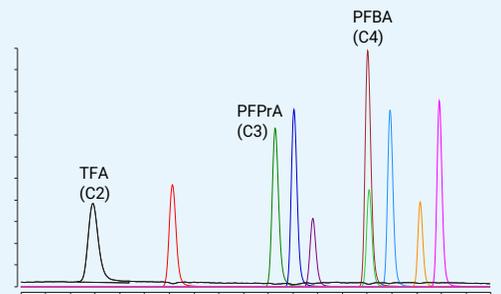
## Engineered for PFAS:

Agilent Altura Poroshell PFAS columns combine a purpose-built stationary phase with advanced Ultra Inert technology—delivering razor-sharp peaks and exceptional consistency, run after run.



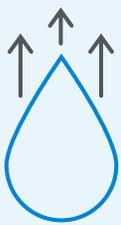
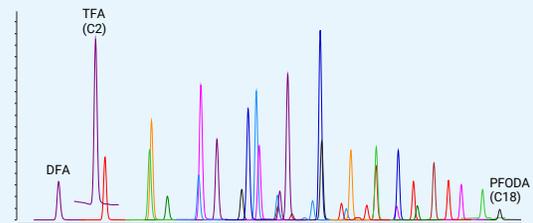
## Boost confidence in every result

- Directly inject drinking water—no sample prep required
- Secure reliable retention for TFA and ultrashort-chain PFAS
- Minimize system background with dedicated delay column



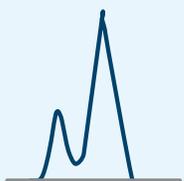
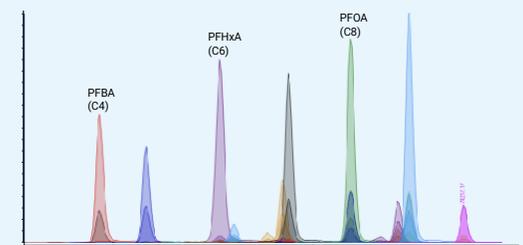
## Accelerate your workflow

- Dilute and shoot C1 to C18 PFAS in a single injection
- Achieve robust, reproducible retention from TFMS and TFA to PFHxDA and PFODA
- Increase throughput with fewer steps, consumables, and reruns



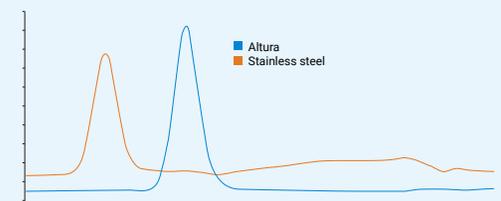
## Inject large sample volumes

- Maintain peak efficiency with high-organic injections up to 200  $\mu$ L
- Ensure consistent retention, even for early eluting PFAS



## Deliver sharper peaks

- Minimize nonspecific binding with Altura Ultra Inert technology
- Support precise, regulatory-ready quantitation across the PFAS panel



Learn more at:  
[www.agilent.com/columns/altura](http://www.agilent.com/columns/altura)

DE.012232

This information is subject to change without notice.

© Agilent Technologies, Inc. 2026  
Published in the USA, January 16, 2026  
5994-8901EN

