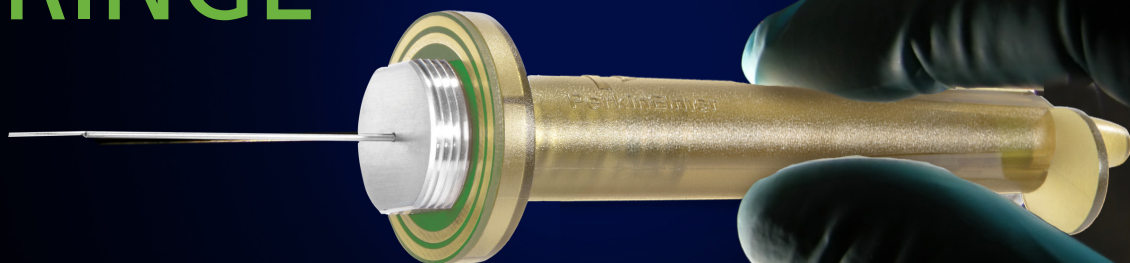


IT'S NOT JUST A SYRINGE



Custodion Sampling Devices

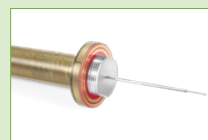
Sample preparation in the field can be difficult with limited resources and time, and your choice of sampling devices can make all the difference. Our rugged Custodion® samplers allow you to achieve quick and reliable results all while on-site with minimal preparation. Our Custodion portfolio includes novel collection and injection devices used for in-field sampling and gas chromatographic (GC) injections. When used with the Torion® T-9 portable GC/MS system, the Custodion devices are a quick, easy, and reliable way to collect and analyze complex samples in the field. Our Custodion line of devices offers an encompassing, flexible solution that meets the diverse and challenging needs of in-field analysis.

The Custodion handles were designed with the field user in mind. They're made of hardened plastic for robust, durable operation, even while wearing heavy protective gear. The push-button trigger on top operates like a ballpoint pen. Each sampler comes with a screw-on/off cap to protect your samples during transport and storage.

The entire Custodion portfolio includes three different sampling devices that provide you with all-encompassing, flexible solutions that meet the diverse, changing needs of sample analysis in the field.

Our range of sampling devices when paired with the Torion T-9 GC/MS offers a complete field-testing package for many different scenarios:

- Rapid identification of chemicals that threaten the public
- Narcotic and drug sample analysis
- Quantitative air quality testing
- Fast screening of water quality



Solid Phase
Microextraction (SPME)



Coiled Microextraction
(CME)

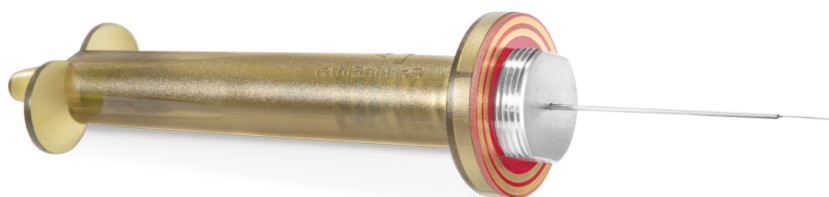


Needle Trap (NT) with
Clairion Pump

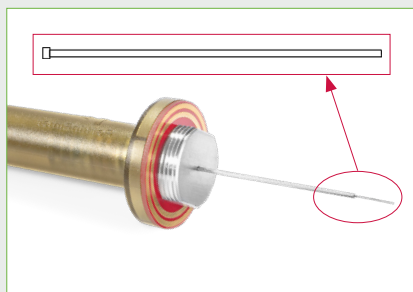


Torion T-9

Custodion Selection Guide



Solid-Phase Microextraction (SPME)



Features:

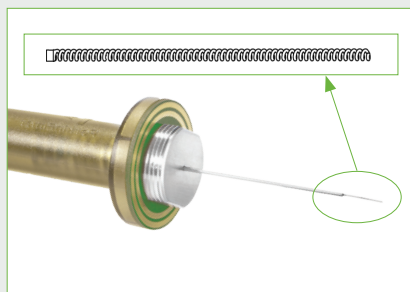
- Passive sampling
- SPME fiber: PDMS / DVB 65 μm
- Concentrates chemical compounds by adsorption
- Headspace or direct fiber immersion

Applications / Ideal Sample Matrix:

- Volatiles, amines and nitro-aromatic compounds (MW 43-300)
- Environmental analysis of water samples
- Forensic analysis of arson/explosive samples
- Flavor analysis of food/beverages

Learn more about our
SPME samplers by visiting
www.perkinelmer.com/...

Coil Microextraction (CME)



Features:

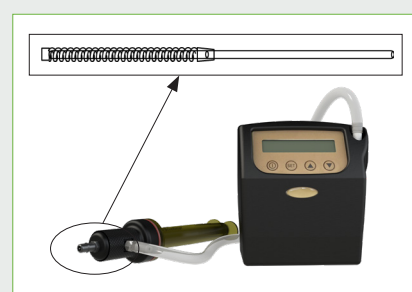
- Inert stainless steel coil draws up dissolved solids and liquid samples
- Chemical contaminants remain on coil when liquid matrix evaporates
- Direct liquid immersion

Applications / Ideal Sample Matrix:

- Semi-volatile, polar compounds (MW 43-500)
- Forensic analysis of pharmaceuticals/drugs of abuse/explosives/CWA
- Essential oil analysis
- Pesticide analysis

Learn more about our
CME samplers by visiting
www.perkinelmer.com/...

Needle Trap (NT)



Features:

- Quantitative analysis with the internal standard module SPS 3
- Active sampling
- Protected tri-bed fiber: Tenax TA, Carboxen 1016 and Carboxen 1003
- Compatible with sample prep station modules: thermal desorption, purge and trap

Applications / Ideal Sample Matrix:

- Trace (ppt) volatile compounds (MW 43-300)
- Environmental and petrochemical analysis of water samples
- Toxic industrial compound analysis
- Air quality monitoring and analysis

Learn more about our
NT samplers by visiting
www.perkinelmer.com/...