Discover the Possibilities

Agilent GC/MS best practices

Perform Sample Cleanup

Sufficient sample preparation and sample cleanup is crucial to getting the most out of any GC/MS analysis.

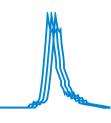


and target

interferences

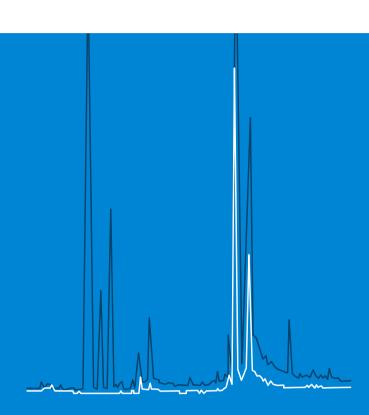


Enhance signal-to-noise ratios and accuracy



Achieve greater reproducibility





Screen Sample Matrix

Performing matrix screening in full scan data acquisition mode facilitates the evaluation of in-source matrix loading.



Prevent source overload



Improve accuracy of analysis



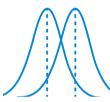
Maintain optimal performance

Dilute More, Inject Less

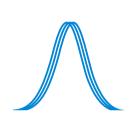
The GC/MS High-Efficiency Source (HES) is designed for ultimate sensitivity where a little goes a long way.



Significantly reduce interferences from matrix background



Peaks with horizontal line



GC/MS operation with excellent ultra-trace level sensitivity





Reduce

Use Backflush

Mid-column backflushing extends the maintenance-free operation of the GC/MS system.



analysis time



Extend column life



Significantly reduce carryover



Quicker inlet and front column maintenance

Implement JetClean

Patented JetClean technology uses a gentle flow of hydrogen through the source.



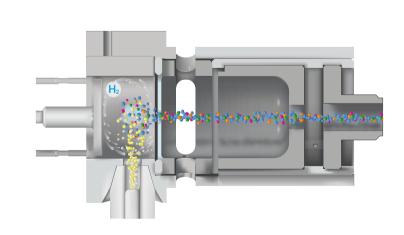
Save time cleaning, reconditioning, and recalibration



Achieve more consistent responses for target analytes



Keep the GC/MS source clean for increased sample throughput



Check for Leaks

To improve your GC/MS analysis, preventing leaks is critical. Agilent offers hardware and software tools to aid in identification of leaks at their source.



The Agilent CrossLab CS Electronic Leak Detector tool detects leaks of various gas types.



Leak test software functionality for GC/TQ identifies the source of leaks in real time when using a leak testing gas.

This information is subject to change without notice.