

Half Century of GC/MS Innovation

Agilent's industry-leading GC/MS systems: The right choice for reproducible results.

1976
5992A Benchtop GC/MS
Up until this time, all GC/MSs were floor-standing units. The 5992 marked a milestone with the first benchtop instrument.



1971
5930A Tabletop MS
The first GC/MS from HP with an oscilloscope and strip chart.



1995 6890 GC 2007 7890A 2016 Intuvo 9000 2019 8860, 8890

1988
Unix and DOS ChemStation
The Unix ChemStation was the successor to the Pascal workstation. The Agilent DOS ChemStation included low-cost PCs and more sophisticated operating systems that made it possible to move to more common computing platforms.



1982
5970 MS
One of Agilent's landmark products, since it was the first of a long line of instruments. With a mass range almost as good as the early floor-standing models, the sensitivity was comparable to our earlier benchtops.



1994
GCD
GC/MS as a technology gained so much popularity that we introduced an easier-to-use model, the GCD.



1996
5973 GC/MSD
The 5973 with positive and negative chemical ionization, showcased extended mass range and sensitivity. MSD ChemStation and the Local Control Panel allowed two GC/MSDs to be controlled by one PC that could be located somewhere else in the lab.



Hyperbolic gold quad
Hyperbolic, gold-coated, quartz quads enhance sensitivity, performance, spectra, and isotope ratios.

2007
MassHunter software
From instrument settings to data analysis and reporting, MassHunter software has made GC/MS analysis both powerful and routine for all.



2009
7000A triple quadrupole GC/MS
The first GC/MS system from Agilent to harness the selectivity and associated sensitivity gains of true MS/MS capability.



2012
7200 GC/Q-TOF
The ideal tool to solve complex problems, the 7200 GC/Q-TOF introduced high resolution accurate mass to Agilent's GC/MS portfolio.



Removable ion source
The removable ion source (RIS) allowed vent-less exchange of EI and CI source technologies on Agilent's 7200 GC/Q-TOF.

2013
5977A GC/MSD
The 5977A introduced the Extractor EI Ion Source for enhanced sensitivity and thermal profile improvement. It also featured direct communication between 7890B GC and the MSD.



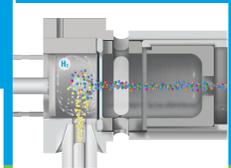
2015
5977B GC/MSD and high efficiency ion source
Delivers unparalleled analytical sensitivity with extreme operational efficiencies for ultra trace-level applications.



2016
7010B triple quadrupole GC/MS
Agilent's 7010B is the evolution of proven performance, featuring compatibility with the high efficiency and JetClean sources, plus the introduction of dMRM Acquisition.



2017
JetClean self-cleaning ion source
JetClean self-cleaning ion source greatly reduces or eliminates the need for source cleaning, thereby enhancing productivity on the single and triple quadrupole GC/MS systems.



7250 GC/Q-TOF
Featuring simultaneous high resolution and high dynamic range, the 7250 enhances and expands upon the hires accurate mass workflows of its predecessor, the 7200 GC/Q-TOF.

2019
QuickProbe GC/MS
The Agilent QuickProbe GC/MS system was designed for forensic drug laboratories facing large sets of samples and growing caseloads.



For more on Agilent GC/MS innovation, visit www.agilent.com/chem/GCMS