

# AGILENT ULTIVO TRIPLE QUADRUPOLE LC/MS

## A SMALL POWERHOUSE OF BIG BENEFITS

### MAXIMIZE LABORATORY REAL ESTATE

The Ultivo LC/TQ makes its mark by reducing the size of its footprint by 70%. Triple your lab's capacity in the same space, with the same power as the big guys.

SMALLER FOOTPRINT

70%

### AGILENT JET STREAM ION SOURCE

More ions created, better MS results

Enhanced desolvation efficiency over a wide range of LC flows.

### VACSHIELD

Optimize lab technician productivity

Enables lab personnel to seamlessly maintain the MS, reducing downtime.

### INTELLIGENT DIAGNOSTICS

Maximum instrument operation to pinpoint issues quickly

Detect issues fast with Ultivo's Intelligent Diagnostics ensuring peak performance.

### CYCLONE ION GUIDE

Maximum performance in a smaller footprint

Get increased signals and more ions to the detector to start producing better results.

### HYPERBOLIC QUADRUPOLE ASSEMBLY

Improved mass resolution and transmission

A truly hyperbolic surface for better mass transmission and improved analytical results.

### VORTEX COLLISION CELL

Increase sample throughput with enhanced MS/MS performance

Efficient collimation of the ion beam with zero crosstalk gives you the confidence to do more faster.

### VIRTUAL PRE/POST-FILTERS

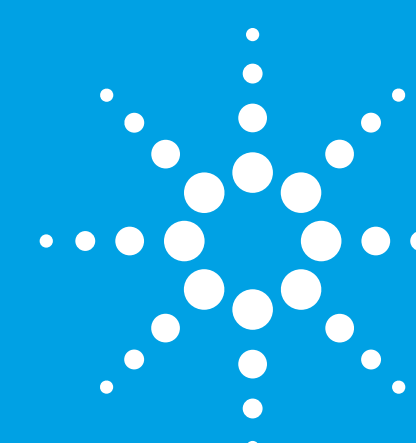
Maximum ion transmission at entrance and exit of resolving quadrupoles

Reduces fringe field effects, allowing for faster switching and enabling more MRMs to be performed.



### ADVANCE EFFICIENCY

The efficiency-optimized Agilent InfinityLab LC series system can sit directly on top of Ultivo, making it the perfect partner for high-throughput mass spectrometry.



Agilent Technologies