



End of Guaranteed Support- LC/MS Systems.

List of LC/MS Systems which reached the End of Guaranteed Support.

Model.	Description.	Introduction.	End of Production.	End-of-Guaranteed-Support.
LC/MSD Single Quad Systems				
<input type="checkbox"/> G1946A	Quadrupole LC/MS	01-Dec-1996	01-Mar-2000	28-Feb -2007
<input type="checkbox"/> G1946B	Quadrupole LC/MS	01-Aug-1999	15-Jun-2000	30-Jun-2007
<input type="checkbox"/> G1946C(VL)	Quadrupole LC/MS	01-Apr-2000	01-Sep-2004	31-Aug-2011
<input type="checkbox"/> G1946D(SL)	Quadrupole LC/MS	01-Apr-2000	01-Sep-2004	31-Aug-2011
<input type="checkbox"/> G1956A	LCMSD Quad VL	01-Jun-2003	01 AUG 2007	30 JUN 2014
<input type="checkbox"/> G1956B	LCMSD Quad SL	01-Jun-2003	01 AUG 2007	30 JUN 2014
<input type="checkbox"/> G6110A	LCMSD Single Quad VL	01-Jun-2006	Current Production	
<input type="checkbox"/> G6120A	LCMSD Single Quad VL Plus	01-Jun-2006	Current Production	
<input type="checkbox"/> G6130A	LCMSD Single Quad SL	01-Jun-2006	Current Production	
<input type="checkbox"/> G6140A	LCMSD High Throughput	01-Jun-2006	Current Production	

This document is believed to be accurate and up-to-date. However, Agilent Technologies, Inc. cannot assume responsibility for the use of this material. The information contained herein is intended for use by informed individuals who can and must determine its fitness for their purpose.

Model.	Description.	Introduction.	End of Production.	End-of-Guaranteed-Support.
LC/MS Trap Systems				
<input type="checkbox"/> G2445A	LCMS Trap	01-Feb-2001	30-Sep-2001	31-Oct-2008
<input type="checkbox"/> G2445C	LCMS Trap VL	01-May-2001	Current Production	
<input type="checkbox"/> G2445D	LCMS Trap SL	01-May-2001	Current Production	
<input type="checkbox"/> G2446A	LCMS Trap XCT	01-May-2003	Current Production	
<input type="checkbox"/> G2447A	LCMS Trap XCT Plus	01-May-2004	Current Production	
<input type="checkbox"/> G2449A	LCMS Trap XCT Ultra	01-May-2005	Current Production	
LC/MSD QQQ Systems				
<input type="checkbox"/> G6410A	LCMSD QQQ	01-Mar-2006	Current Production	
<input type="checkbox"/> G6410B	LCMSD QQQ	01-Mar-2008	Current Production	
LC/MSD QTOF Systems				
<input type="checkbox"/> G6510A	LCMSD QTOF	01-Aug-2003	Current Production	
<input type="checkbox"/> G6520A	LCMSD QTOF	01-Sep-2007	Current Production	
LC/MSD TOF Systems				
<input type="checkbox"/> G1969A	LCMSD TOF	01-Aug-2003	Current Production	
<input type="checkbox"/> G6220A	LCMSD TOF	01-Sep-2007	Current Production	

This document is believed to be accurate and up-to-date. However, Agilent Technologies, Inc. cannot assume responsibility for the use of this material. The information contained herein is intended for use by informed individuals who can and must determine its fitness for their purpose.

www.chem.agilent.com

This information is subject to change without notice.

© Agilent Technologies, Inc. 2007
Printed in USA August 7, 2007