



Agilent Technologies

Ordering Agilent 1200 Series Rapid Resolution LC System

The Agilent 1200 Series Rapid Resolution LC (RRLC) System can be ordered in various configurations to meet the different application and performance requirements. Three configuration examples are described below, which span the full performance and functionality range of the Agilent 1200 RRLC system.

Whatever system configuration is chosen as a starting point - the modular, scalable and open system architecture of the Agilent 1200 RRLC system allows future expandability and adaptability to other configurations and applications. Please contact your local sales representative for more detailed information and further advice.

Agilent 1200 Series RRLC System - Configuration Examples

Configuration 1

A cost-effective configuration offering full compatibility with Rapid Resolution LC, Ultra-fast LC and conventional HPLC on standard bore columns with 3 - 4.6 mm id, 10 - 300 mm length and particle sizes from 1.5 to 10 μm . This configuration includes the Agilent 1200 Series VWD SL for high-speed single-wavelength UV-Vis detection

Configuration 2

The most flexible standard configuration offering full compatibility with Rapid Resolution, Ultra-fast and conventional LC and LC/MS on narrow and standard bore columns with 2.1 - 4.6 mm id, 10 - 300 mm length and particle sizes from 1.5 to 10 μm . This configuration includes the DAD SL for highest-speed multi-wavelength and full spectral UV-Vis detection for peak purity analysis and spectral conformation.

Configuration 3

A sophisticated high-throughput configuration with alternating column regeneration (ACR) that offers ultimate throughput of up to 2000 samples per day - in a sequential manner without loss in performance or method flexibility. This configuration provides full compatibility with Rapid Resolution, Ultra-fast and conventional LC and LC/MS on narrow and standard bore columns with 2.1 - 4.6 mm id, 10 - 300 mm length and particle sizes from 1.5 to 10 μm . It includes the Agilent 1200 Series DAD SL for highest-speed multi-wavelength and full spectral UV-Vis detection for peak purity analysis and spectral conformation.

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.

Agilent 1200 Series RRLC System - Configuration 1

Description	Product Number	Price and Availability
Agilent 1200 Series micro vacuum degasser	G1379B	Not available online
Agilent 1200 Series binary pump SL	G1312B	Not available online
Agilent 1200 Series high performance autosampler SL	G1367C	Not available online
Agilent 1200 Series thermostatted column compartment SL	G1316B	Not available online
Agilent 1200 Series variable wavelength detector SL	G1314C	Not available online
Standard flow cell for Agilent 1200 Series VWD SL 10-mm pathlength, 14- μ L volume, pressure tight to 40 bar	G1314C #018	Not available online
Semi-micro flow cell for Agilent 1200 Series VWD SL 6-mm pathlength, 5- μ L volume, pressure tight to 40 bar	G1314C #016	Not available online

Agilent 1200 Series RRLC System - Configuration 2

Description	Product Number	Price and Availability
Agilent 1200 Series micro vacuum degasser	G1379B	Not available online
Agilent 1200 Series binary pump SL	G1312B	Not available online
Seal wash option for Agilent 1200 Series binary pump SL	G1312B #030	Not available online
Solvent selection option for Agilent 1200 Series binary pump SL	G1312B #031	Not available online
Agilent 1200 Series high performance autosampler SL	G1367C	Not available online
Agilent 1200 Series thermostat for high performance autosampler SL	G1330B	Not available online
Agilent 1200 Series thermostatted column compartment SL	G1316B	Not available online
Low dispersion kit Capillaries and heat exchangers for compatibility with 2.1mm narrow bore columns	G1316B#060	Not available online
Agilent 1200 Series diode-array detector SL	G1315C	Not available online

Standard flow cell for Agilent 1200 Series diode-array detector SL 10-mm pathlength, 13- μ L volume, pressure tight to 120 bar	G1315C #018	Not available online
Semi-micro flow cell for Agilent 1200 Series diode-array detector SL 6-mm pathlength, 5- μ L volume, pressure tight to 120 bar	G1315C #016	Not available online
Micro flow cell for Agilent 1200 Series diode-array detector SL 3-mm pathlength, 2- μ L volume, pressure tight to 120 bar	G1315C #010	Not available online
Agilent 1200 Series RRLC System - Configuration 3		
Description	Product Number	Price and Availability
Agilent 1200 Series micro vacuum degasser	G1379B	Not available online
Agilent 1200 Series binary pump SL	G1312B	Not available online
Seal wash option for Agilent 1200 Series binary pump SL	G1312B #030	Not available online
Solvent selection option for Agilent 1200 Series binary pump SL	G1312B #031	Not available online
Agilent 1200 Series Binary Pump SL Second Pump for alternating column regeneration	G1312B	Not available online
Seal wash option for Agilent 1200 Series Binary Pump SL	G1312B #030	Not available online
Agilent 1200 Series high performance autosampler SL	G1367C	Not available online
Agilent 1200 Series thermostat for high performance autosampler SL	G1330B	Not available online
Agilent 1200 Series thermostatted column compartment SL	G1316B	Not available online
2-position/10-port micro valve SL for Agilent 1200 Series thermostatted column compartment SL For alternating column regeneration. Includes valve and accessory kit.	G1316B#057	Not available online
Low dispersion kit Capillaries and heat exchangers for compatibility with 2.1mm narrow bore columns	G1316B#060	Not available online
Alternating column regeneration (ACR) kit Capillaries for alternating column regeneration	G1316B#061	Not available online

Agilent 1200 Series diode-array detector SL	G1315C	Not available online
Standard flow cell for Agilent 1200 Series diode-array detector SL 10-mm pathlength, 13- μ L volume, pressure tight to 120 bar	G1315C #018	Not available online
Semi-micro flow cell for Agilent 1200 Series diode-array detector SL 6-mm pathlength, 5- μ L volume, pressure tight to 120 bar	G1315C #016	Not available online
Micro flow cell for Agilent 1200 Series diode-array detector SL 3-mm pathlength, 2- μ L volume, pressure tight to 120 bar	G1315C #010	Not available online