

# Agilent 1120 Compact LC

## Application Notes

Part Number	Title
5989-7455EN	Analysis of herbicides in drinking water using the Agilent 1120 Compact LC
5989-7456EN	Analysis of antioxidants in chewing gum using the Agilent 1120 Compact LC
5989-7457EN	Development and validation of a method for simultaneous determination of paracetamol, diclofenac and ibuprofen using the Agilent 1120 Compact LC
5989-7458EN	Analysis of ginseng and American ginseng using the Agilent 1120 Compact LC
5989-8330EN	In-process control for Aripiprazole manufacturing with the Agilent 1120 Compact LC and ZORBAX C-18 column
5989-8330ENA	In-process control for Aripiprazole manufacturing with the Agilent 1120 Compact LC and Agilent TC C18(2) columns
5989-8331EN	Detector linearity testing for Aripiprazole quality control with the Agilent 1120 Compact LC and ZORBAX C-18 columns
5989-8331ENA	Detector linearity testing for Aripiprazole quality control with the Agilent 1120 Compact LC and Agilent TC C18(2) columns
5989-8333EN	System suitability testing for Aripiprazole quality control with the Agilent 1120 Compact LC and ZORBAX C-18 columns
5989-8333ENA	System suitability testing for Aripiprazole quality control with the Agilent 1120 Compact LC and Agilent TC C18(2) columns
5989-8334EN	System suitability testing for Tramadol quality control with the Agilent 1120 Compact LC and ZORBAX C-18 columns
5989-8334ENA	System suitability testing for Tramadol quality control with the Agilent 1120 Compact LC and Agilent TC C18(2) columns
5989-8335EN	Injector linearity testing for Tramadol quality control with the Agilent 1120 Compact LC and ZORBAX C-18 columns
5989-8335ENA	Injector linearity testing for Tramadol quality control with the Agilent 1120 Compact LC and Agilent TC C18(2) columns
5989-8960EN	Analysis of preservatives in food and cosmetics with the Agilent 1120 Compact LC
5989-8961EN	Analysis of sweeteners in food and beverages with the Agilent 1120 Compact LC
5989-9241EN	Development and validation of an HPLC method to analyze ibuprofen and impurities according to the European Pharmacopoeia using the Agilent 1120 Compact LC
5989-9869EN	USP purity analysis of pravastatin sodium using the Agilent 1120 Compact LC
5989-9949EN	Quantitative Liquid Chromatography Analysis of Melamine in Dairy Products Using Agilent's 1120 Compact LC and 1200 Rapid Resolution LC