



## CONFERENCE AGENDA

### Automation in the Chromatography Laboratory

Tuesday 6<sup>th</sup> & Wednesday 7<sup>th</sup> October 2009

Down Hall Country House Hotel

Day 1 – Tuesday 6<sup>th</sup> October 2009

11.30 – 12.50	Registration and Lunch	
12.50 – 13.00	Welcome	<i>Mr. Ray Perkins, Anatune Ltd, UK</i>
13.00 – 13.40	The Successful Application of Stir Bar Sorptive Extraction (Twisted) for the Trace Measurement for Different Classes of Compounds in Water	<i>M. David Benanou, Veolia Water, France</i>
13.40 – 14.20	New Technology for Air Analysis and Soil Gas Measurements	<i>Dr. Daniel Cardin, Entech Instruments Inc, USA</i>
14.20 – 15.00	Automated Sample Preparation Using the MPS Autosampler with MAESTRO Software Control	<i>Mr. Fredrick Foster, GERSTEL Inc, USA</i>
15.00 – 15.30	Tea/Coffee Break	
15.30 – 16.10	Considering Automated Sample Preparation in the Context of LC-QQQ	<i>Mr. John Lee, Agilent Technologies, UK</i>
16.10 – 16.50	Automating Dispersive Solid Phase Clean-Up (QuEChERS) for LC-MS and GC-MS	<i>Mr. Paul Roberts Anatune Ltd, UK</i>
18.45 – 19.30	Pre-Dinner Drinks in the Priory Room	
19.30	Dinner	

Day 2 – Wednesday 7<sup>th</sup> October 2009

07.00 – 08.45	Breakfast (served between these times)	
09.20 – 10.00	Practical Applications of Disposable Pipette Extraction (DPX) in Forensic Toxicology	<i>Dr. Simon Elliot, ROAR Forensics, UK</i>
10.00 – 10.40	Molecularly Imprinted Polymers - Highly Selective SPE for Trace Analysis in Complex Matrices	<i>Dr. Lisa Fitzpatrick, Sigma Aldrich, UK</i>
10.40 – 11.20	Hyphenating Automated Solid Phase Extraction with GC-MS	<i>Mr. Jonathan Angove, Anatune Ltd, UK</i>
11.20 – 11.40	Tea/Coffee Break	
11.40 – 12.20	Automated Sample Preparation for LC-MS in the Clinical Laboratory	<i>Dr. Meike Baden, TeLA, Germany</i>
12.20 – 13.00	Automated Sample Preparation in Food Analysis	<i>Dr. Oliver Lerch, GERSTEL GmbH, Germany</i>
13.00 - 13.40	The Use of Stir Bar Sorptive Extraction Back Extraction for the Clean-Up of Clinical Samples Prior to LC-MS	<i>Dr. Prashant Kole The Queen's University, Belfast</i>
13.40	Lunch	

