

Ion Chromatography Workshop

Altrincham, UK | May 17, 2017

Agenda

08:50	Registration
09:20	Welcome and introduction
09:30	Suppression Unravelled: Cost Saving, Sensitivity and Robustness
10:30	Column Considerations: Improve Resolution, Speed of Analysis and Longevity
11:30	Coffee break
11:45	Novel Screening Method Using Dual Channel Ion Chromatography <i>Alastair Bewsher, The University of Manchester</i>
12:15	Sample Preparation: Protect your Consumables, Instrument and Achieve Better Analysis
13:00	Lunch
13:45	5 Quick Tips to Improve or Develop Your Ideal Method
14:15	Carbohydrate Analysis using HPAE-PAD (ICS4000) for Evaluating the Inhibition of Sugar Digestion in an Intestinal Cell Model <i>Alison Pyner, The University of Leeds</i>
14:45	Chromeleon: Tips and Tricks from Start-up to Final Report
15:15	Coffee break
15:30	ULS: A Simple Guide to First Line Troubleshooting
16:00	IC The Future: The Intelligent Laboratory
16:30	Meeting close

Venue:

Thermo Fisher Scientific
1 St George's Court
Altrincham
Cheshire
WA14 5TP