



LC-MS Annual Meeting

London 28 June 2017

Agenda

Registration and coffee 09:00		
09:30	Welcome and Introduction <i>Phil Green, Thermo Fisher Scientific</i>	
09:40	ASMS 2017 Review - Routine Applications <i>Claudia Martins, Thermo Fisher Scientific</i>	
10.05	Precision Mass Spectrometry – ASMS Update <i>Ken Miller, Thermo Fisher Scientific</i>	
10:30	Plenary Lecture Mass Spectrometry of Membrane Proteins and their Complexes <i>Professor Dame Carol Robinson, University of Oxford</i>	
Coffee break 11:15		
Omics Parallel Session		Small Molecules Parallel Session
One Year with a Thermo Scientific™ Orbitrap Fusion™ Tribrid™ Mass Spectrometer <i>Adam Dowle, York University</i>	11:45	Multi-residue analysis of pesticides in water by solid phase extraction coupled to liquid chromatography Orbitrap Mass Spectrometry <i>Jorge Casado, Greenpeace Science Unit</i>
Novel insight to the causal link between insulin signaling deregulation & its potential etiology of aggressive prostate cancer: Implications to the development of novel biomarkers & treatment targets <i>Spiros Garbis, University of Southampton</i>	12:05	Thermo Scientific™ Q Exactive™ for Met ID in a DMPK Lab <i>Lloyd King, UCB</i>
Lunch 12:25		

Venue:

[Wellcome Collection](#)

183 Euston Road

London

NW1 2BE

Afternoon Sessions		
Omics Parallel Session		Small Molecules Parallel Session
An in vitro MS assay for sensitive determination of donor specific MHC class II immunopeptidomes <i>Martin Pabst, Abzena</i>	13:40	Title tbc <i>Allan Traynor, Alere</i>
Glycoproteomics and Intact MS analysis of HIV immunogens <i>Weston Struwe, Oxford Glycobiology</i>	14:00	Novel approaches for toxicology and explosives detection <i>Melanie Bailey, University of Surrey</i>
Lipidomics supporting predictive toxicology – finding the camel in a sandstorm <i>Aniko Kende, Syngenta</i>	14:20	Small Molecule Software - ASMS Update <i>Mark Sanders, Thermo Fisher Scientific</i>
Coffee break 14:40		
Omics Parallel Session		Small Molecules Parallel Session
Get to know your mAb <i>David Muñoz, Thermo Fisher Scientific</i>	15:10	Identification of non-intentionally added substances during migration analysis of printing inks <i>Marina Santos, Domino</i>
Systems biology of oncogenic signalling <i>Pedro Cutillas, Barts Cancer Institute</i>	15:30	The Analysis of Emerging Polar Contaminants by Hyphenated IC-MS/MS <i>Michal Godula, Thermo Fisher Scientific</i>
Large Molecule Software - ASMS Update <i>Mark Sanders, Thermo Fisher Scientific</i>	15:50	New achievements and perspectives in high-performance mass spectrometry imaging <i>Bernhard Spengler, TransMIT GmbH</i>
Imaging Lipid Remodelling in Cancer and Metabolic Disease <i>Zoe Hall, University of Cambridge</i>	16:10	The combined use of the Thermo Scientific™ TSQ Quantiva™ and Q Exactive for enhanced DMPK Discovery support <i>Iain Beattie, XenoGesis</i>
Close and wine reception 16:30		

Find out more at thermofisher.com/uk-ireland-events