

# Tecan Symposium comes to Boston



The fifth annual Tecan Symposium will be visiting the historical US coastal city of Boston, Massachusetts, from the 23<sup>rd</sup> to the 25<sup>th</sup> of October. Building on the success of previous events, this popular scientific meeting is being hosted by Tecan in the US for the first time, and will bring together key opinion

leaders from around the world to discuss the expanding role of mass spectrometry (MS) in life sciences and diagnostics.

This year's scientific program will cover a broad range of topics related to MS, and will be divided into four sessions: the role of MS in basic research; collection, preparation, and storage of samples for MS; MS as a diagnostic tool; and the use of MS in applied markets. Supported by a number of social events, this diverse agenda is designed to attract a truly multidisciplinary audience. As in previous years, places at the Symposium are limited

to ensure plenty of opportunities for open discussions and networking, giving speakers and delegates the chance to explore novel concepts, applications and solutions that might help to address unmet scientific needs, while enjoying Tecan's hospitality.

To find out more about the Tecan Symposium, including details regarding free registration, visit [www.tecan.com/symposium](http://www.tecan.com/symposium)

## Tecan Symposium key speakers

### Prof Sabine Becker

BrainMet: novel avenues of bioimaging mass spectrometry of metals and biomolecules from micrometer to nanometer scale in the brain

### Prof Mitsutoshi Setou

Meta-analysis of histopathological lipidomics

### Dr Feixia Chu

Sample preparation of ancient protein samples from amber encapsulated organisms

### Dr Gary van Berkel

Laser ablation and liquid extraction surface sampling for ambient surface sampling/ionization MS

### Dr Jeffrey Hurst

Chocolate spectrometry: applications of mass spectrometry in cocoa and chocolate research

### Dr Michael Vogeser

Pitfalls associated with the use of liquid chromatography-tandem mass spectrometry in the clinical laboratory

### Prof Xuefan Gu

Neonatal screening using tandem mass spectrometry

### Dr John Brennan

Mass spectrometric methods for biological screening of small molecule mixtures

### Dr Nigel Clarke

Mass spectrometry as an enabling technology in the clinical laboratory

### Dr Graham Bench

Emerging trends and technologies enabling ultrasensitive <sup>14</sup>C measurements for biomedicine

### Dr Paul Tempst

Aminopeptidase activities as biomarkers for cancer

### Dr Judith Stone

Automating high throughput LC-MSMS analysis of 25-hydroxy vitamin D using liquid handlers and middleware

### Dr Robert Moritz

The human SRMAtlas: a compendium of quantitative mass spectrometry assays for all human proteins

### Dr Andrew Hoofnagle

Developing methods for LC-MS-based clinical diagnostics

### Dr Mark Libardoni

The role of mass spectrometry in solar system geochemistry and planetary atmospheres: past missions to state-of-the-art instrument development

### Dr Burak Eral

eMALDI: suppressing the coffee stain effect via electrowetting for improved MALDI-MS detection

### Dr Zoltan Takats

Mass spectrometric profiling of biological tissues: a comprehensive alternative to classical histopathology

### Dr Stuart Black

The application of stable isotope radio mass spectrometry for forensic analysis of human skeletal remains

### Dr Mehdi Moini

Biological clocks: high throughput identification of deterioration markers and dating of museums' proteinaceous specimens