

Structural mass spectrometry for the analysis of complex biological samples: ion mobility-mass spectrometry for broad-scale systems and synthetic biology

John A. McLean

**Laboratory for Structural Mass Spectrometry
Department of Chemistry
Vanderbilt Institute for Chemical Biology
Vanderbilt Institute for Integrative Biosystems Research and Education
Vanderbilt University
Nashville, TN 37235**

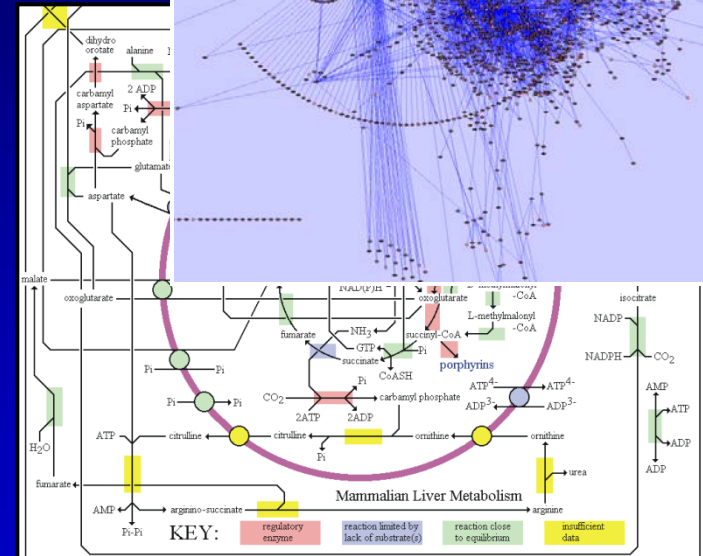
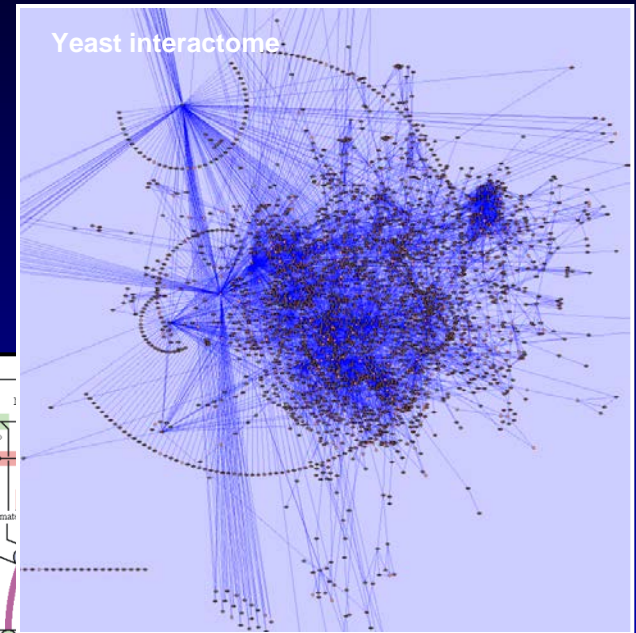


**GEN – Genetic Engineering and Biotechnology News
March 25th, 2014**

Integrating omics for systems biology

Characterizing, quantifying, and cataloging the biomolecular inventory of a sample at specific dimensions of:

- (i) Space (e.g. cellular, tissue, or organism level)
- (ii) Time (e.g. point in the life cycle, healthy vs. diseased state).



One of the grand challenges in Systems Biology is the development of new technologies for characterizing the dynamic temporal change of the biomolecular inventory...

Paraphrasing Leroy Hood
Pittcon Heritage Award Address 2008

On the dimensions of cartography...

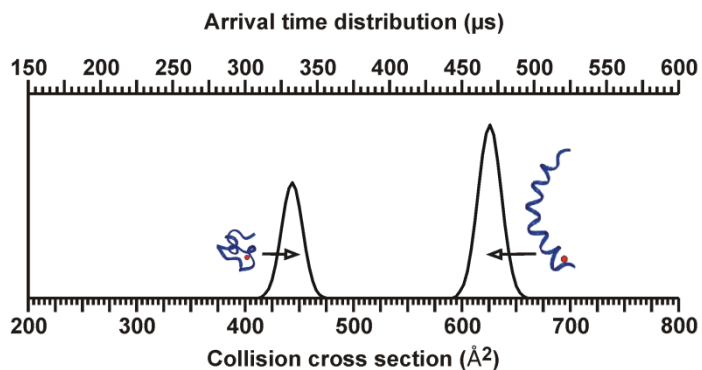
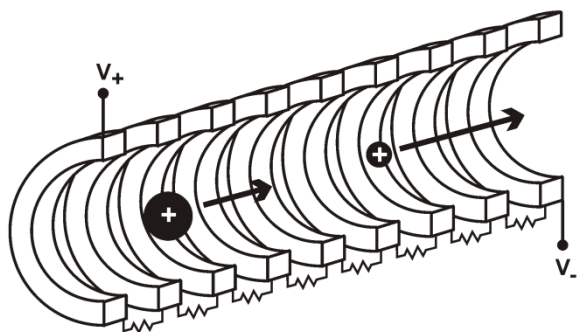
Suarez Miranda, Viajes de varones prudentes, Libro IV, Cap. XLV, Lerida, 1658:

“. . . In that Empire, the Art of Cartography attained such Perfection that the map of a single Province occupied the entirety of a City, and the map of the Empire, the entirety of a Province. In time, those Unconscionable Maps no longer satisfied, and ***the Cartographers Guilds struck a Map of the Empire whose size was that of the Empire, and which coincided point for point with it.*** The following Generations, who were not so fond of the Study of Cartography as their Forebears had been, saw that that vast Map was Useless, and not without some Pitilessness was it, that they delivered it up to the Inclemencies of Sun and Winters. In the Deserts of the West, still today, there are Tattered Ruins of that Map, inhabited by Animals and Beggars; in all the Land there is no other Relic of the Disciplines of Geography.”

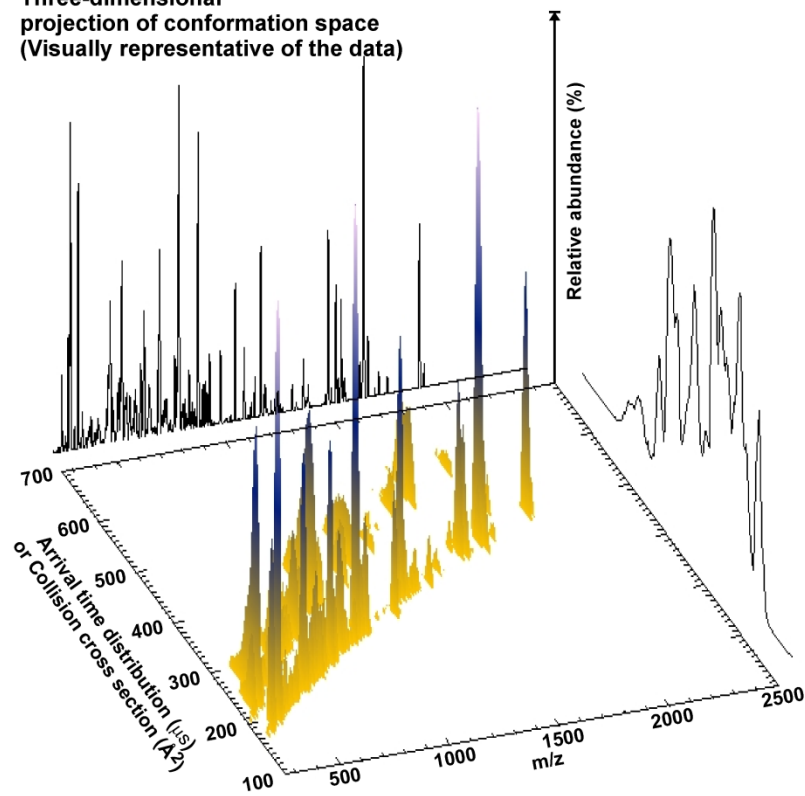
(From Jorge Luis Borges, Collected Fictions translated by Andrew Hurley)

On the dimensions of dynamics in omics...
A broad timescale ranging from genomics to metabolomics

On the dimensions of ion mobility-mass spectrometry...



Three-dimensional projection of conformation space (Visually representative of the data)



Molecular structure determines collision cross section

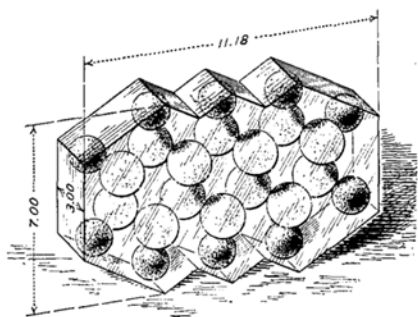
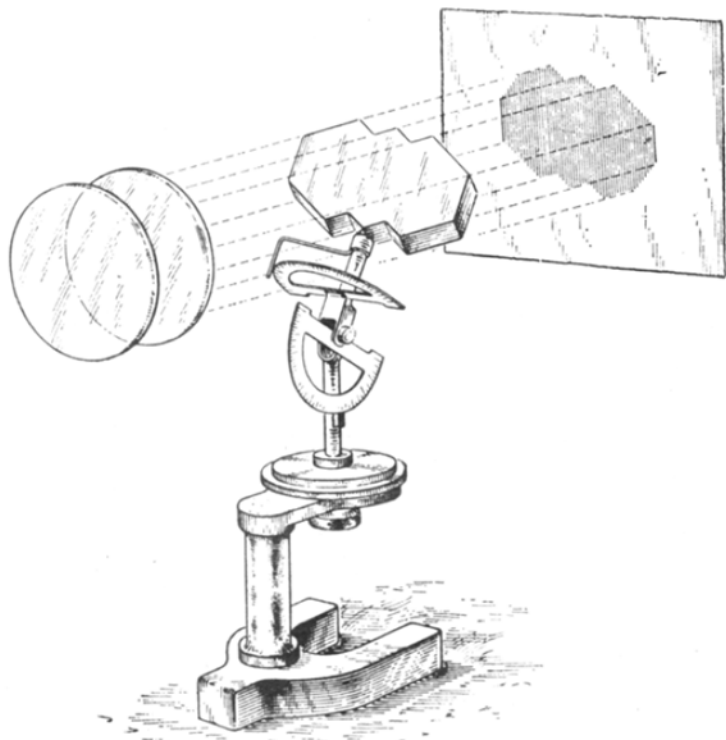


Fig. 7.—Anthracene.

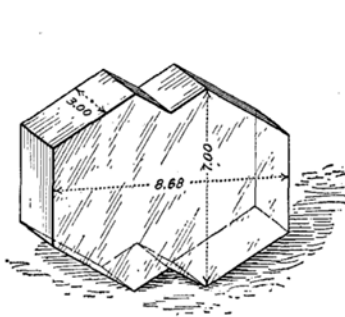
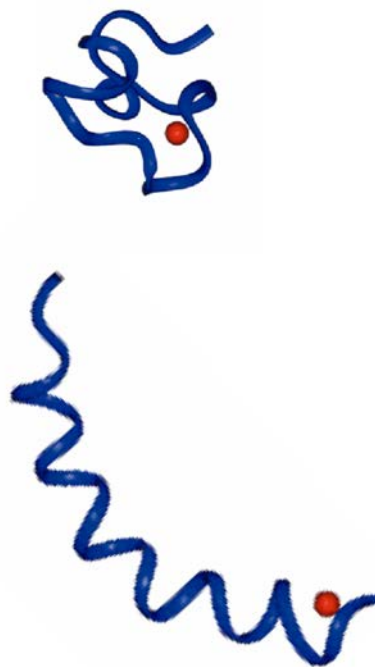
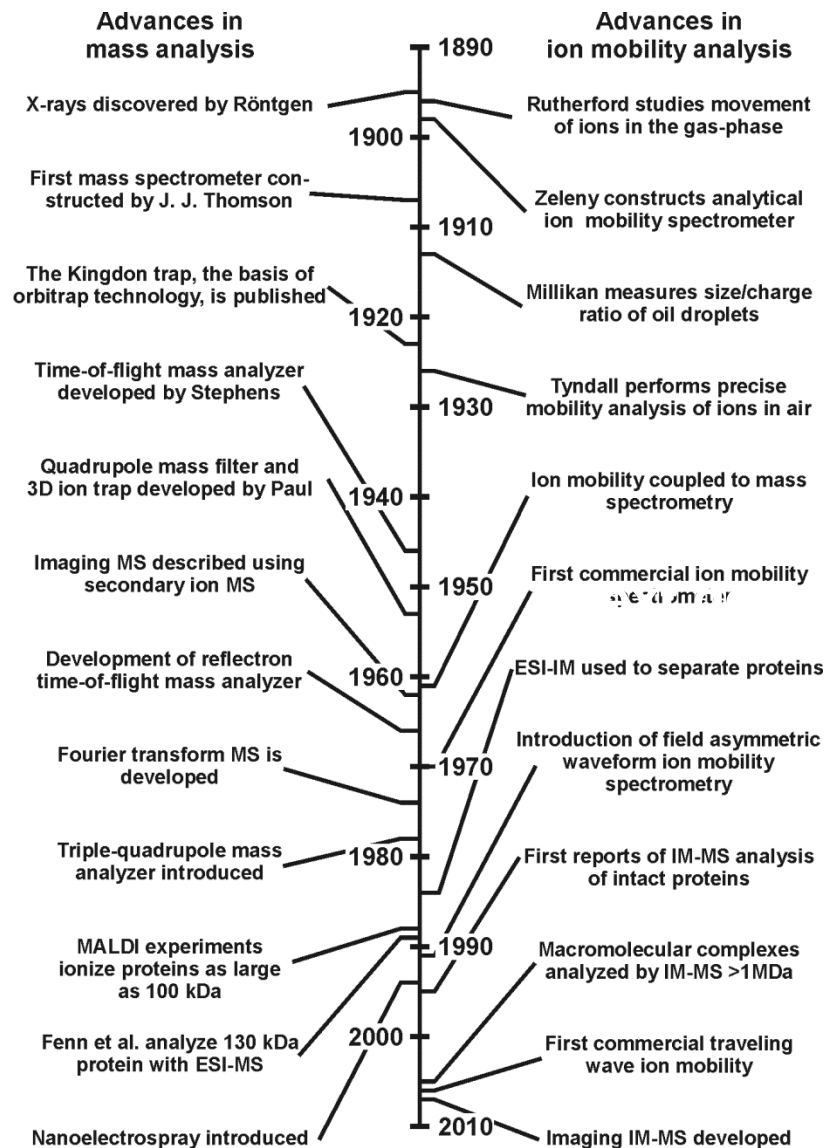


Fig. 8.—Naphthalene.



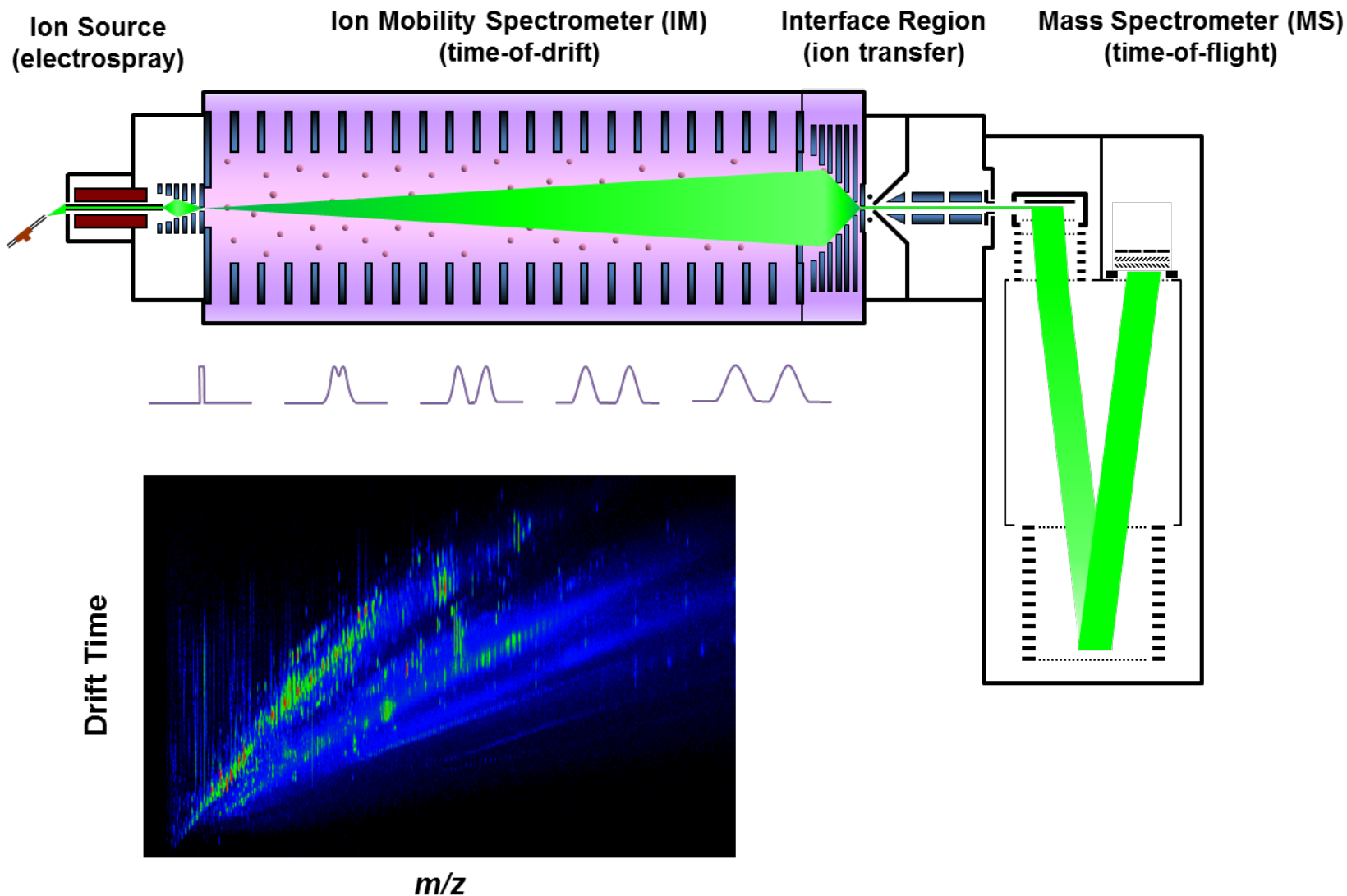
E. Mack Jr., *J. Am. Chem. Soc.* 47, 2468-82 (1925).
Ref. 1. Langmuir, *J. Am. Chem. Soc.* 39, 1848 (1917).
Ref. 3. Bragg, *Proc. Phys. Soc.* 34, 33 (1921).

Evolution of ion mobility with mass spectrometry



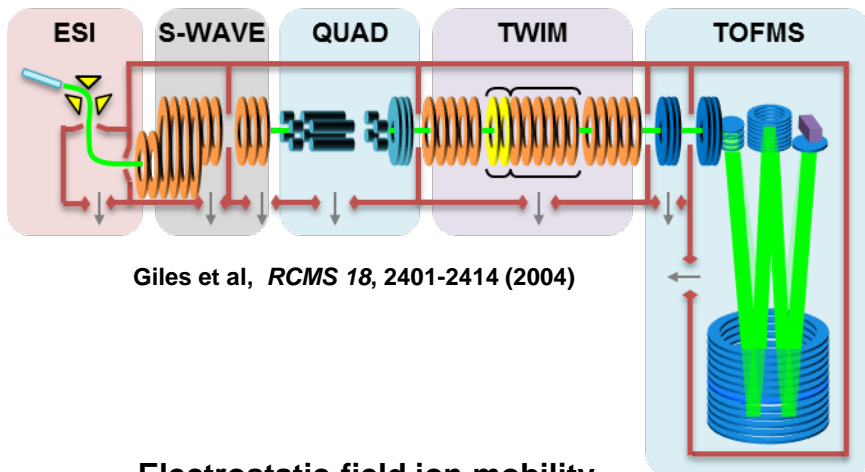
Time-dispersive ion mobility-mass spectrometry

Charge-to-size and mass-to-charge separations



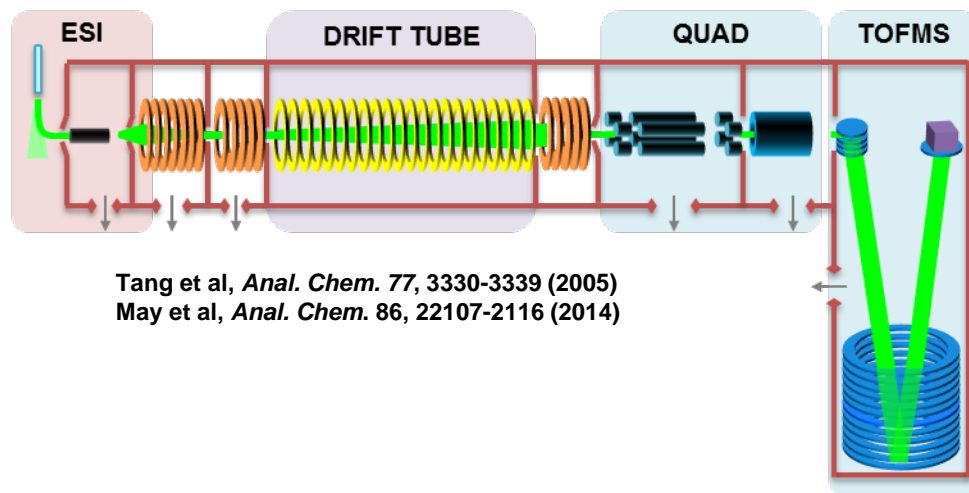
Three major types of time-dispersive ion mobility-mass spectrometry instrumentation

Electrodynamic-field ion mobility
(relative collision cross section)



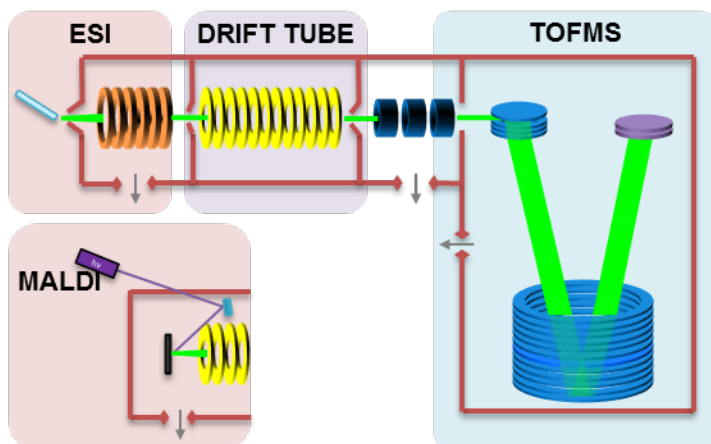
Giles et al, *RCMS* 18, 2401-2414 (2004)

Electrostatic-field ion mobility
(absolute collision cross section)

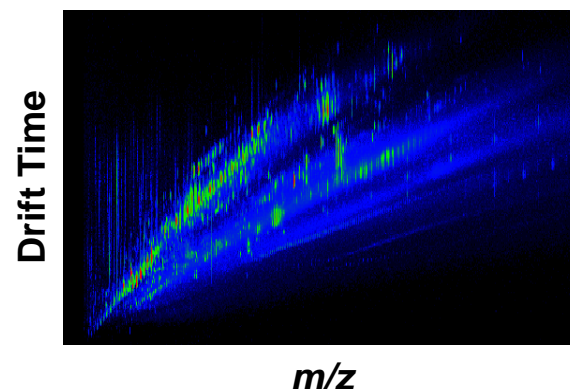


Tang et al, *Anal. Chem.* 77, 3330-3339 (2005)
May et al, *Anal. Chem.* 86, 22107-2116 (2014)

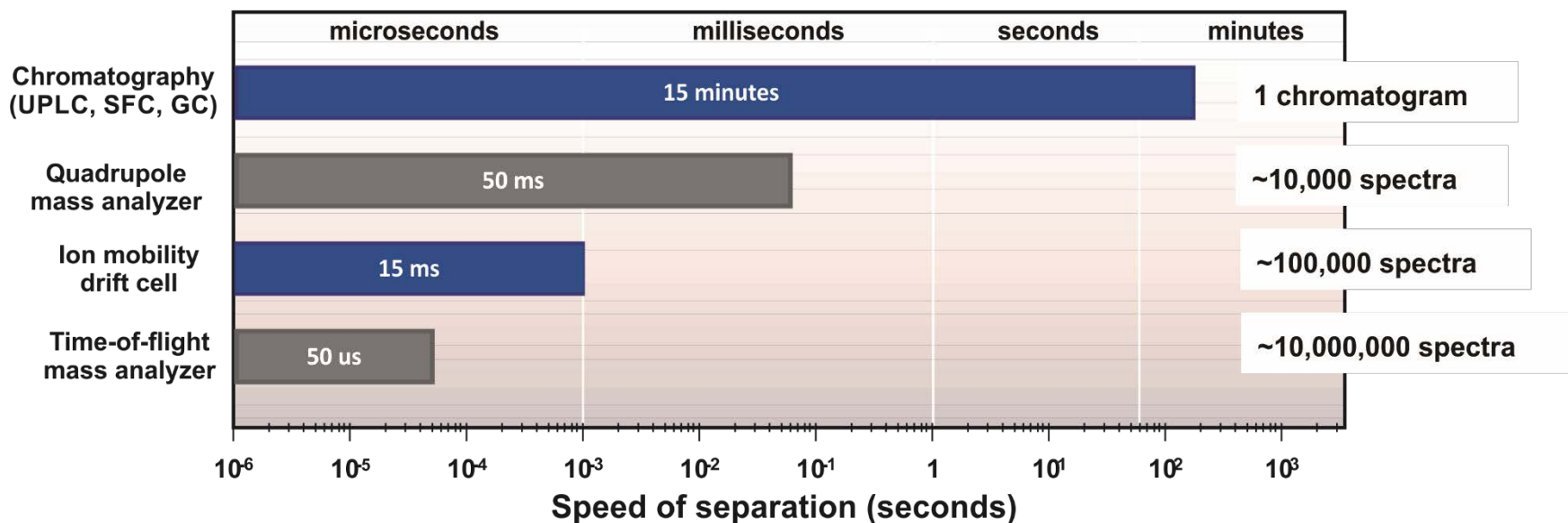
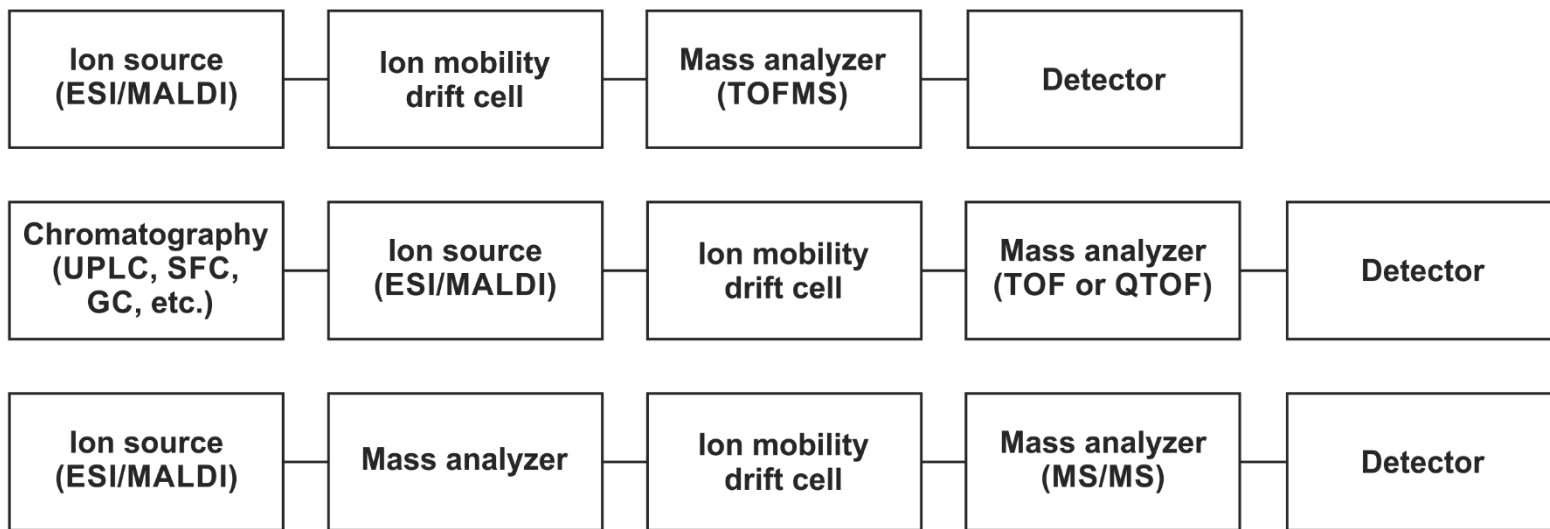
Electrostatic-field ion mobility
(absolute collision cross section)



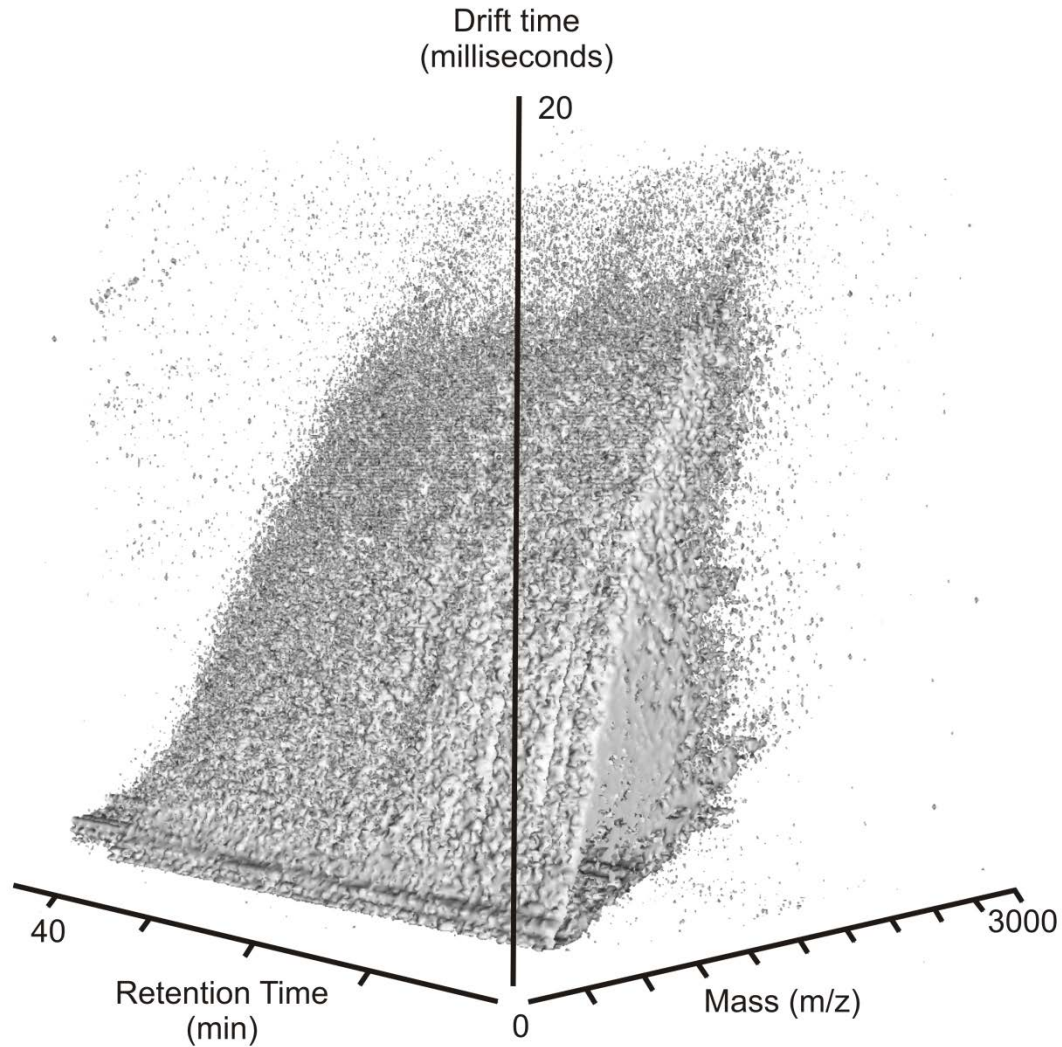
Sundarapandian et al, *Anal. Chem.* 82, 3247-3254 (2010)



Nesting of analytical timescales and data scaling

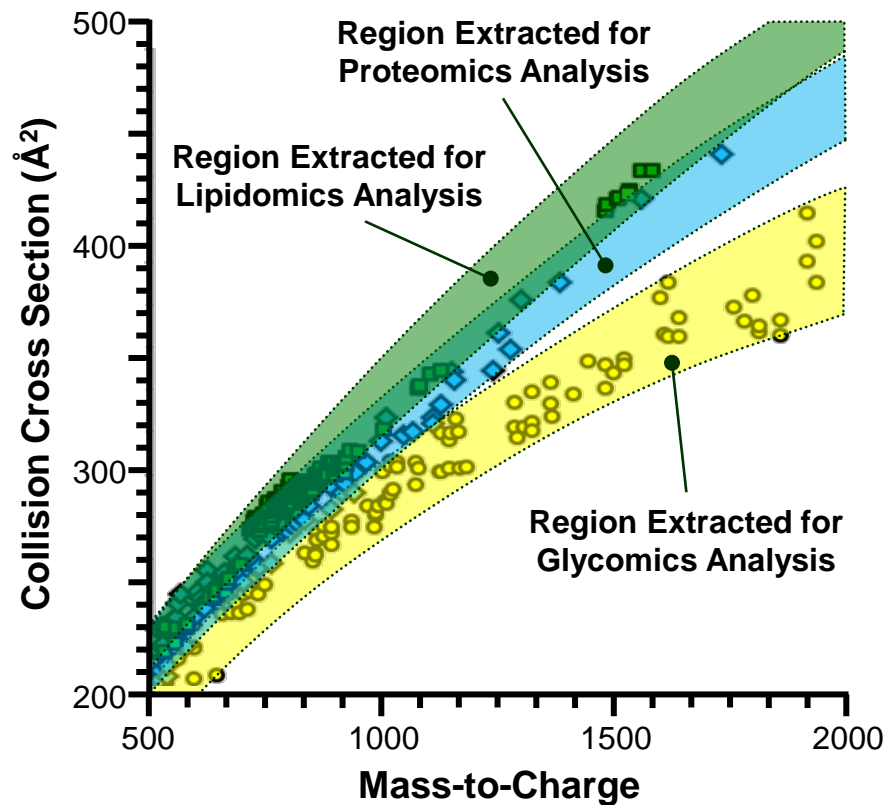
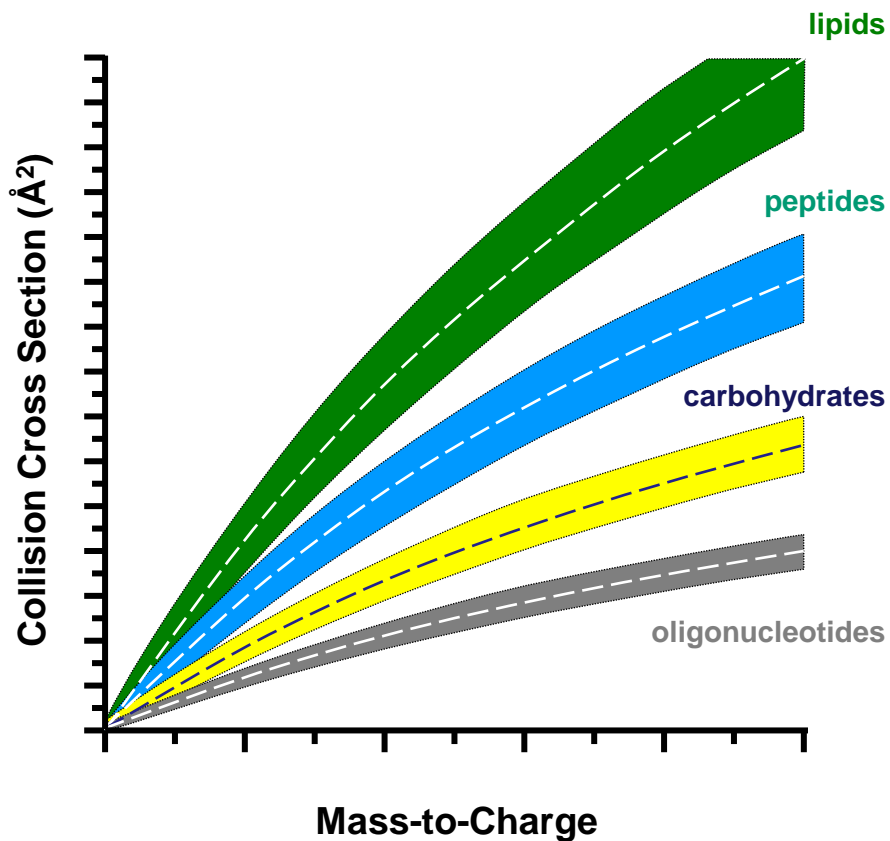


UPLC-IM-MS analysis of serum



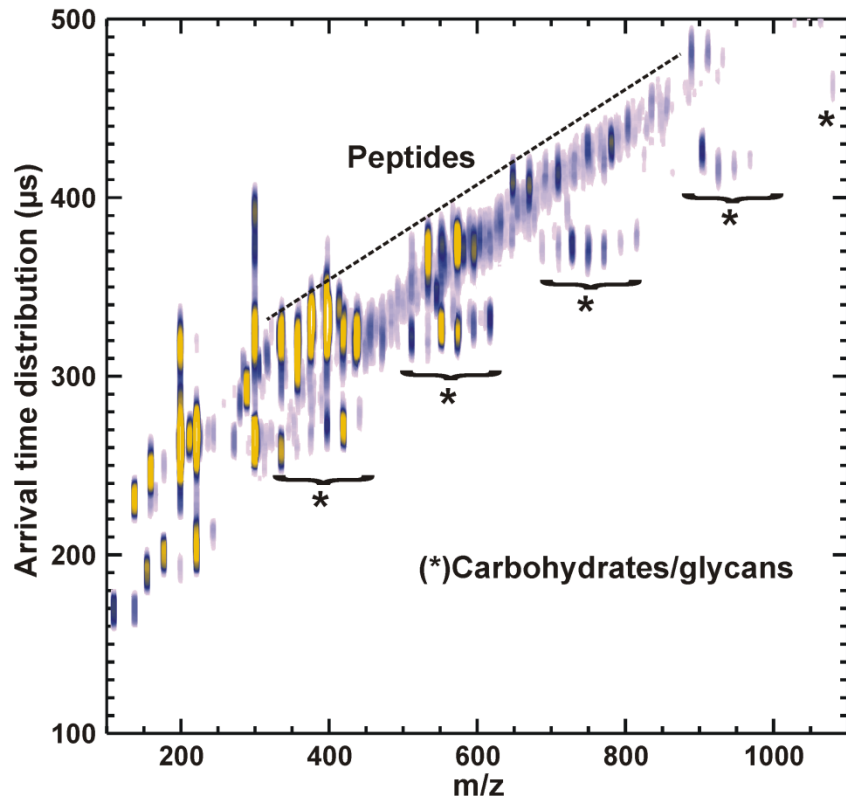
Biomolecular class separations

(Structural selectivity from prevailing structural characteristics of density)

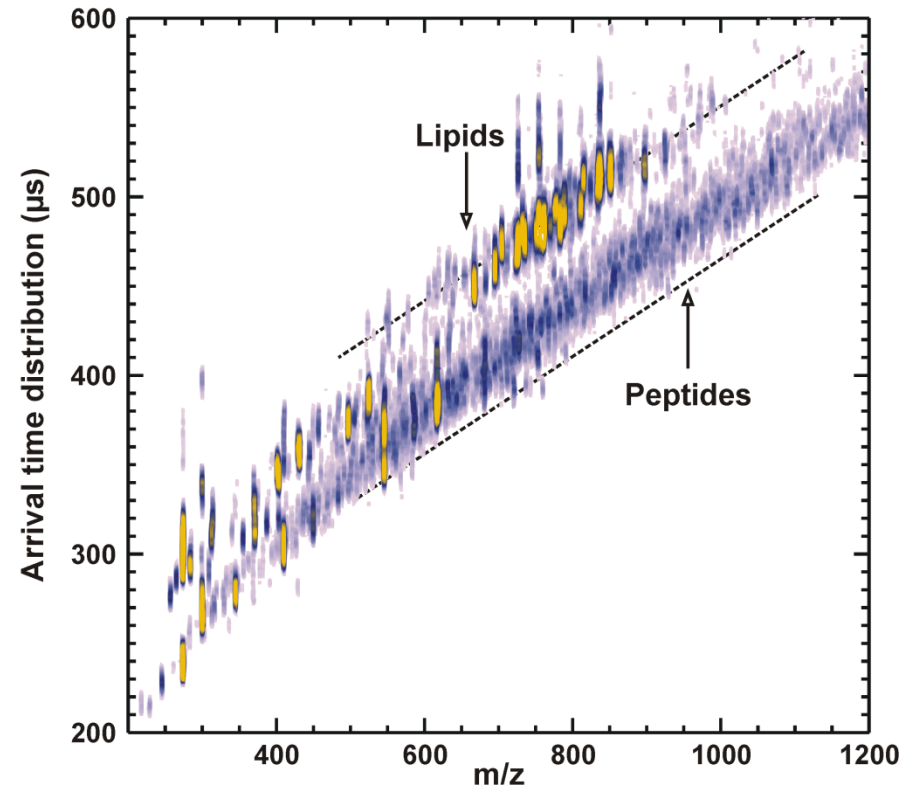


Integrated omics

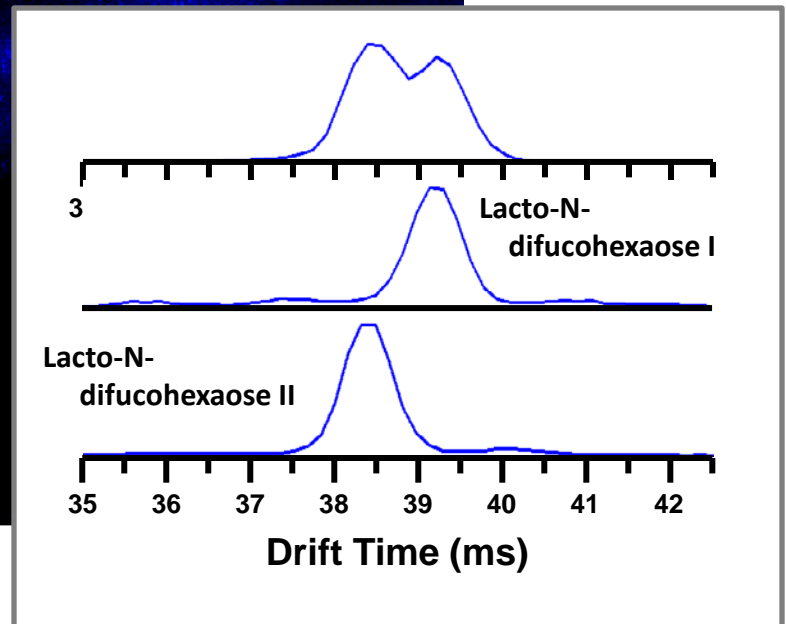
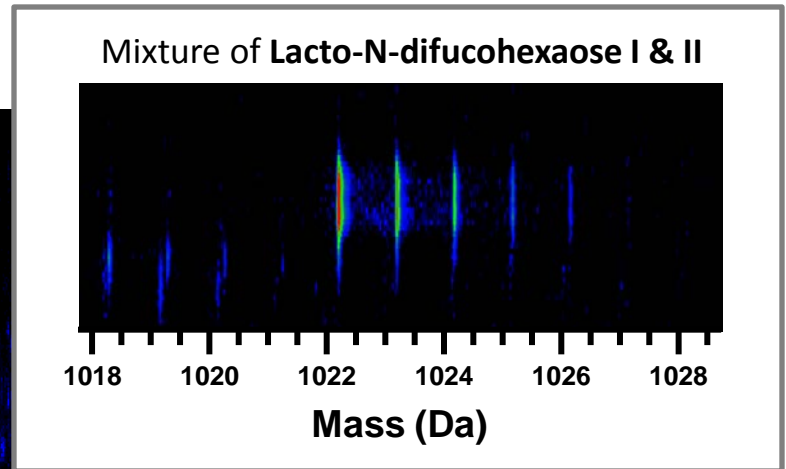
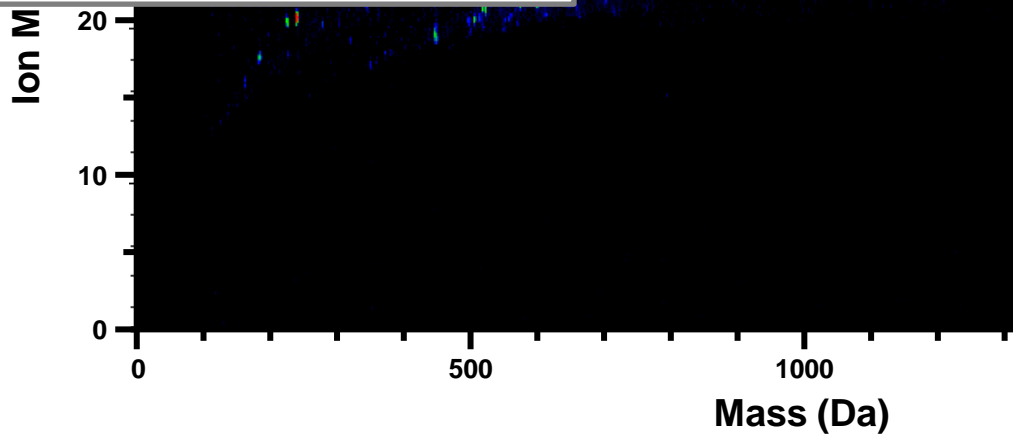
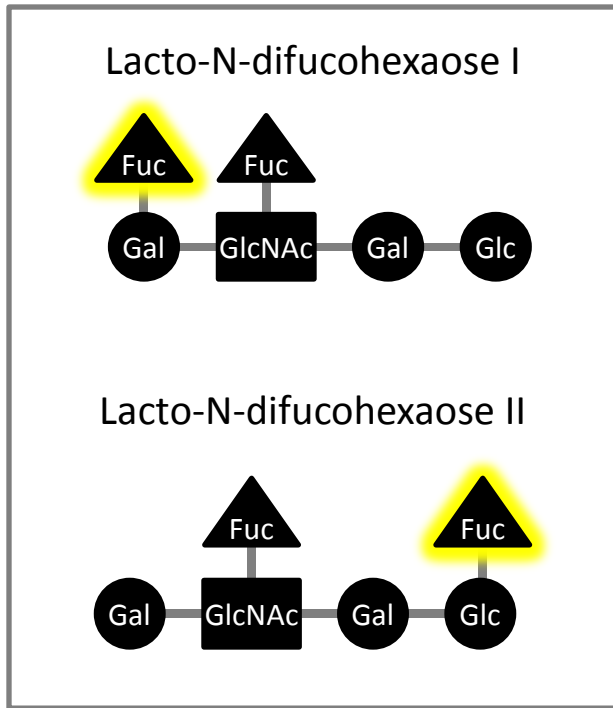
**Simultaneous glycoproteomics
(glycans and peptides from RNase b)**



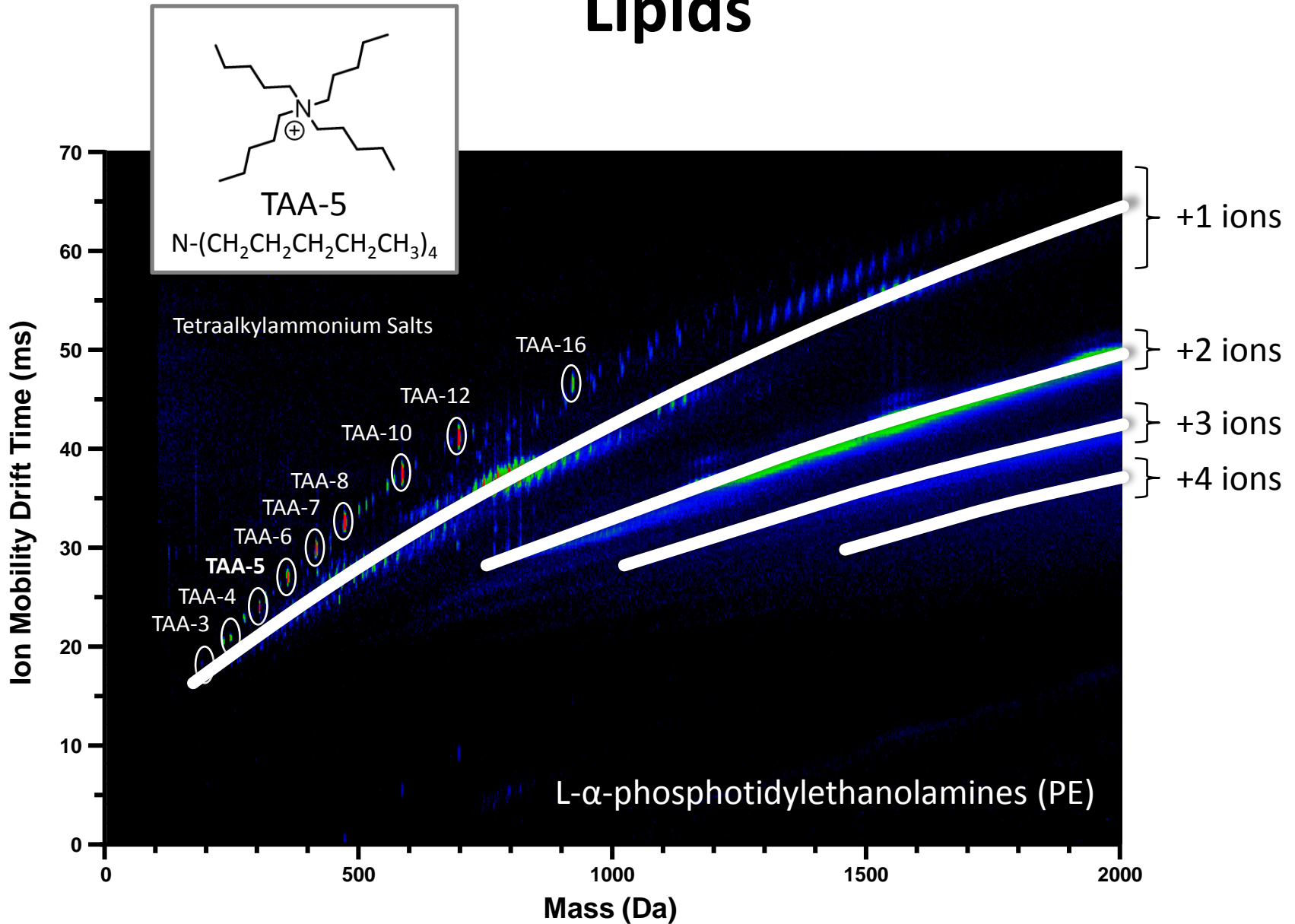
**Simultaneous peptide and lipid profiling
(human glioblastoma)**



Carbohydrates

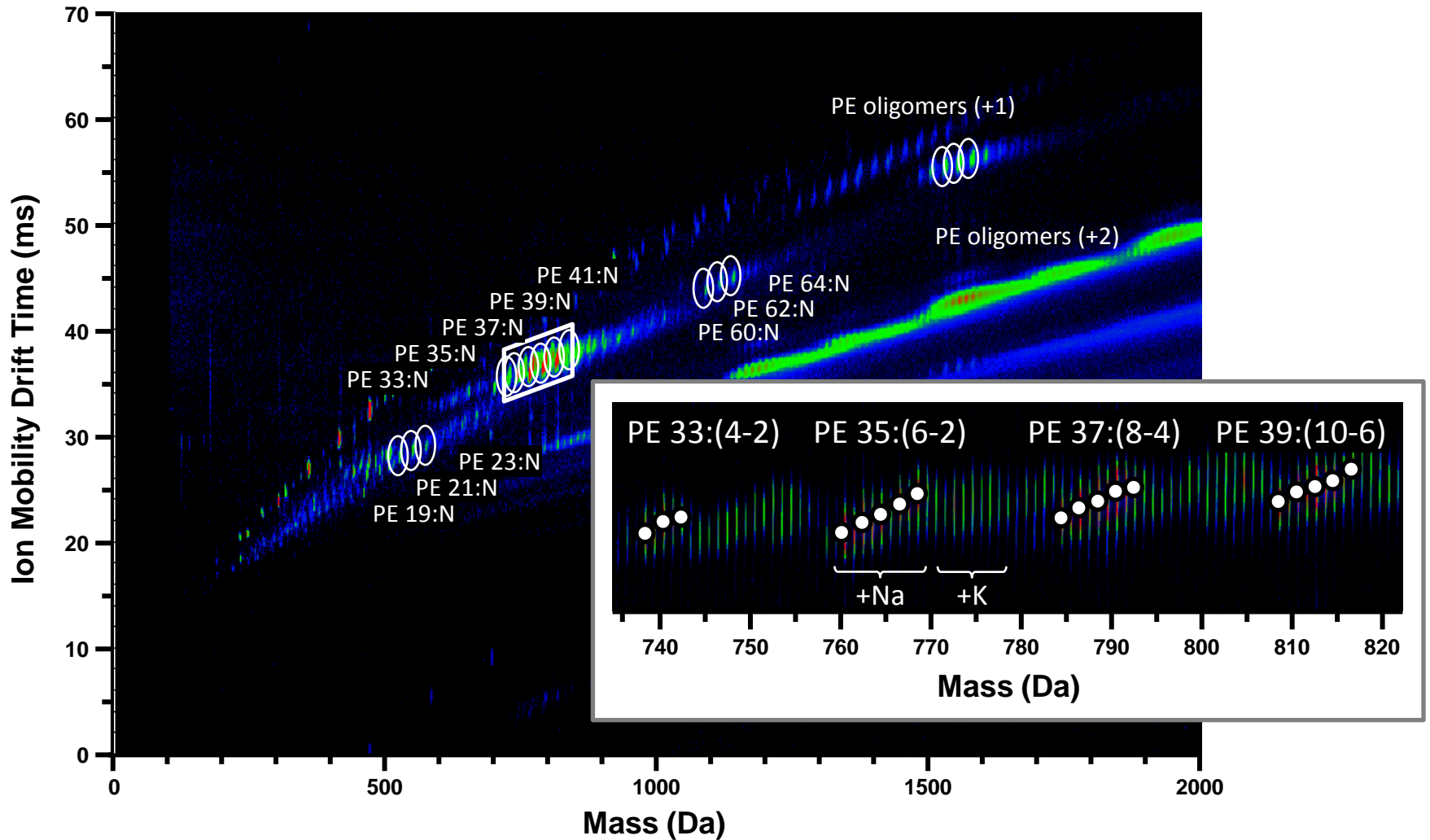


Lipids



Distribution of PE lipids derived from chicken egg. An equimolar mixture of TAA salts was spiked in as an internal IM and MS calibrant. Structure shown is the tetrapentylammonium cation (TAA-5).

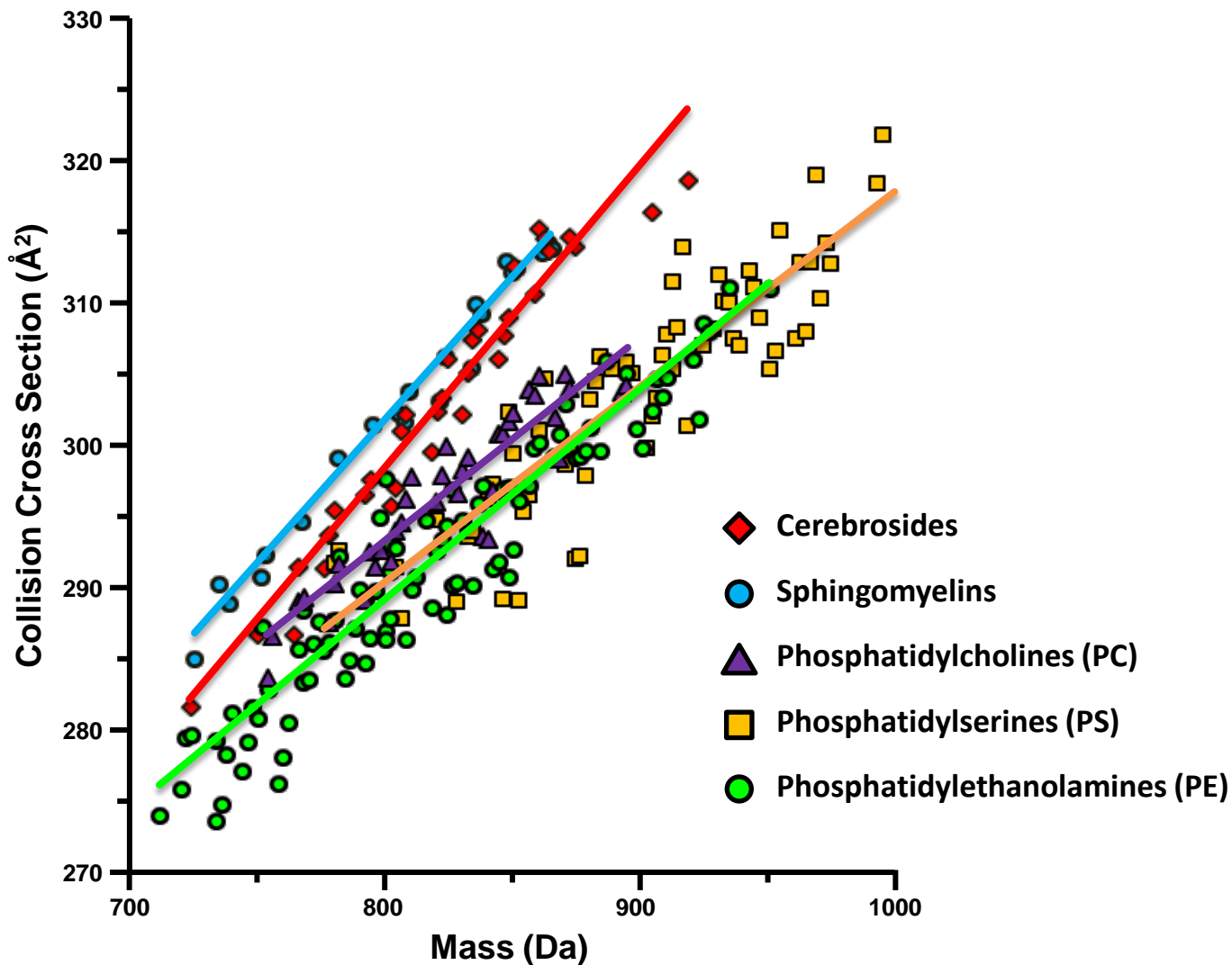
Lipids



Distribution of PE lipids derived from chicken egg. An equimolar mixture of TAA salts was spiked in as an internal IM and MS calibrant.

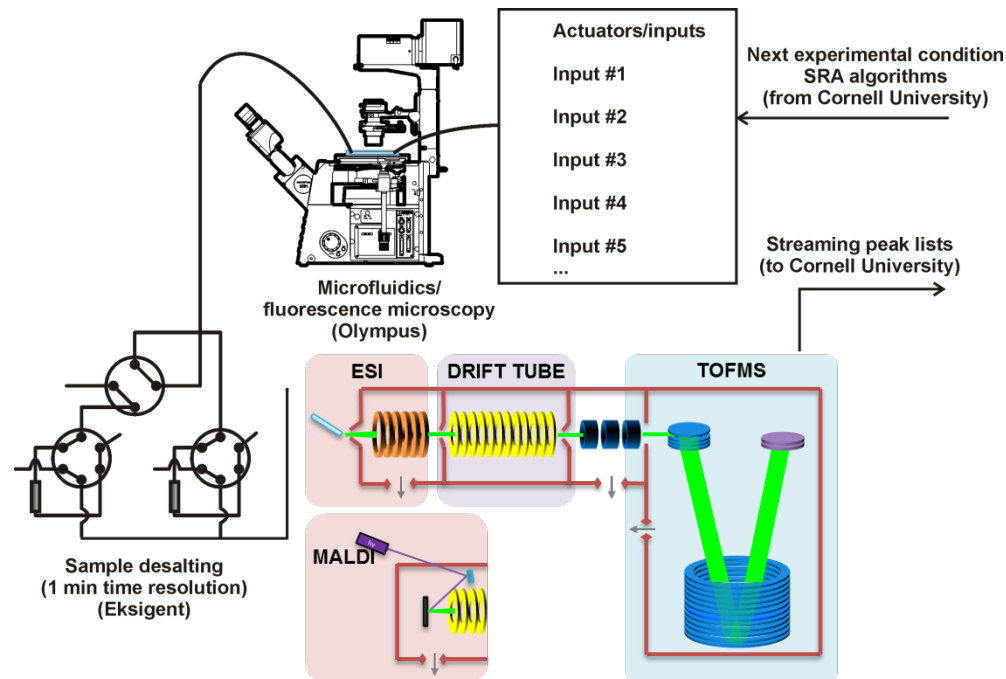
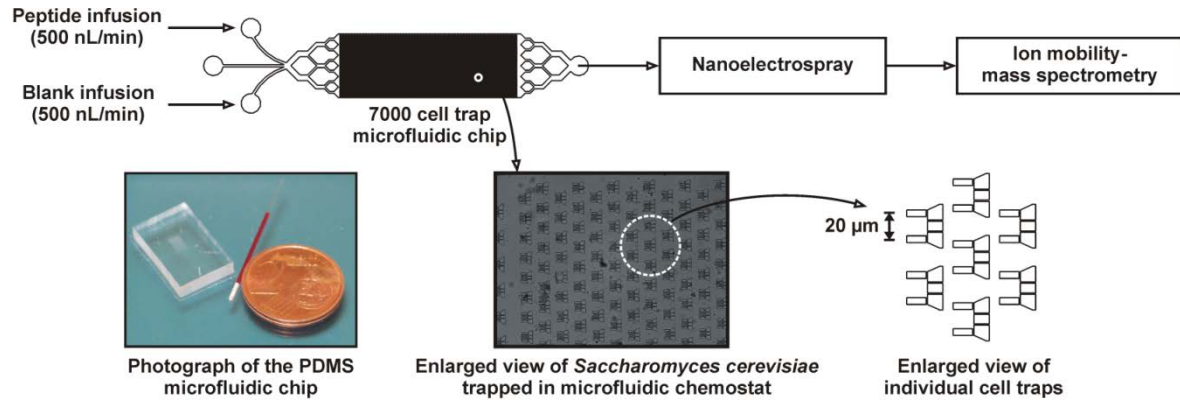
Conformational Space Mapping in High Resolution

Sub-Class Trends in Lipid Data



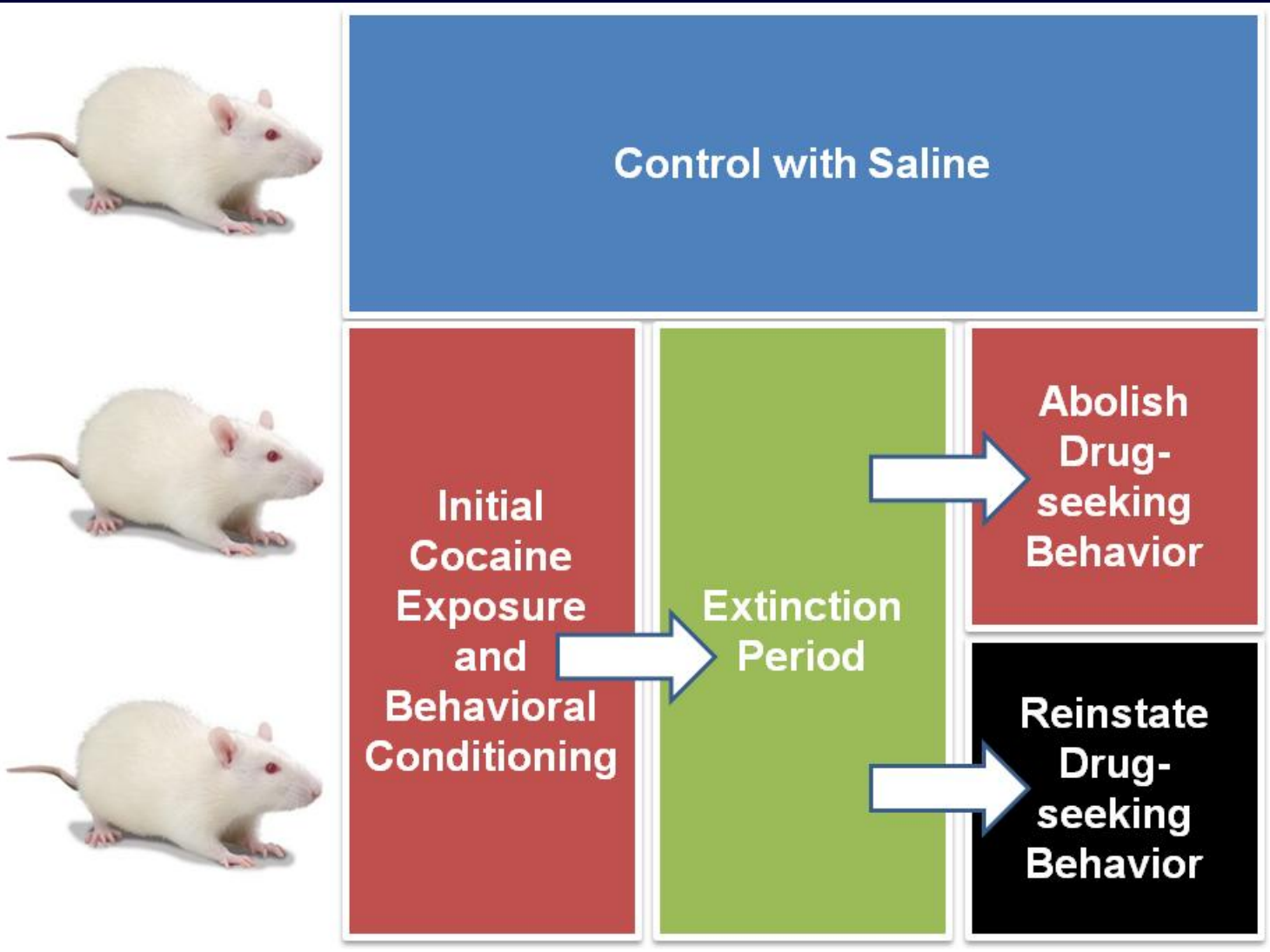
Dynamic systems biology

Automated robot scientist version 2.0

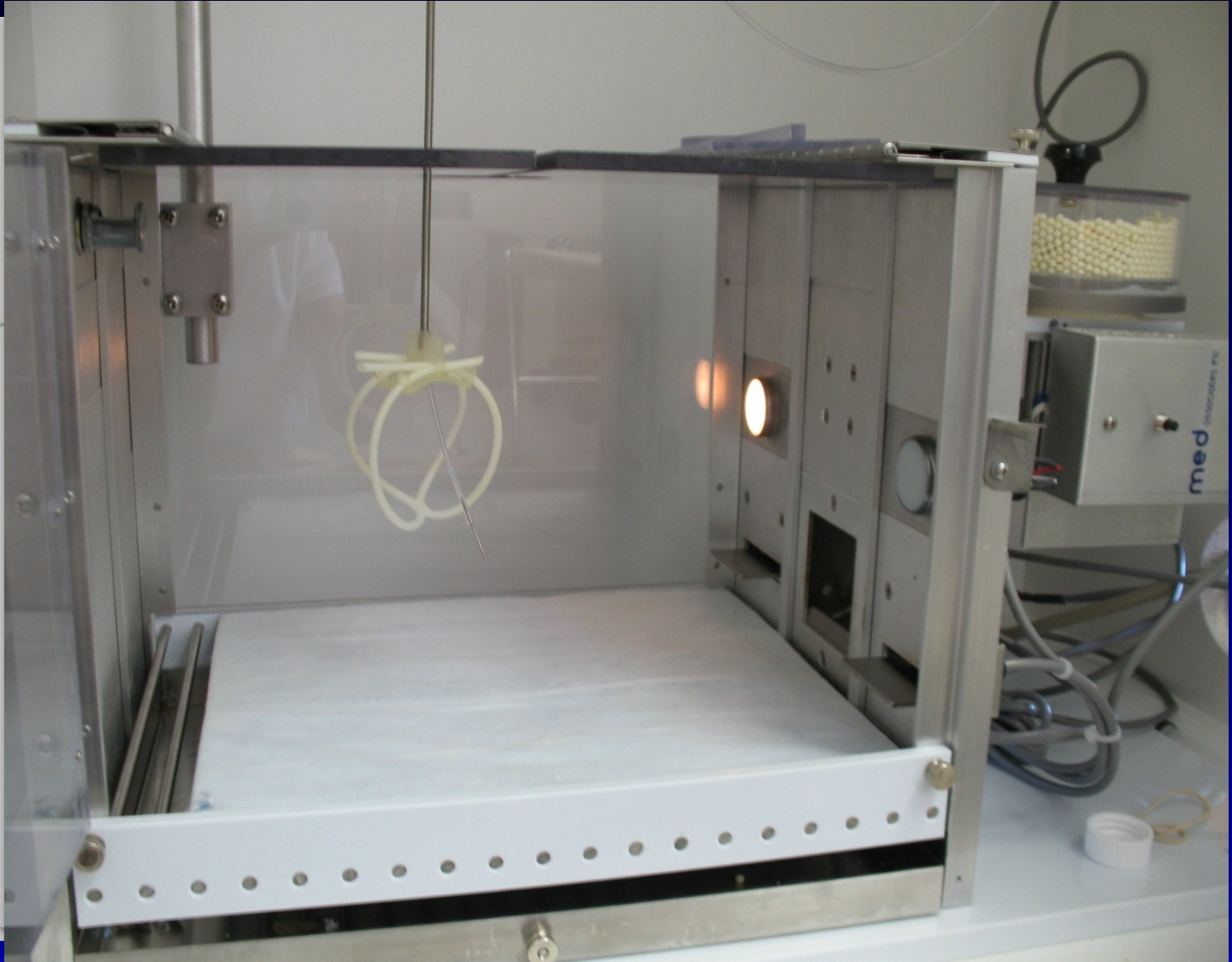


Collaboration with
Prof. John Wikswo (VU) and
Prof. Hod Lipson (Cornell)

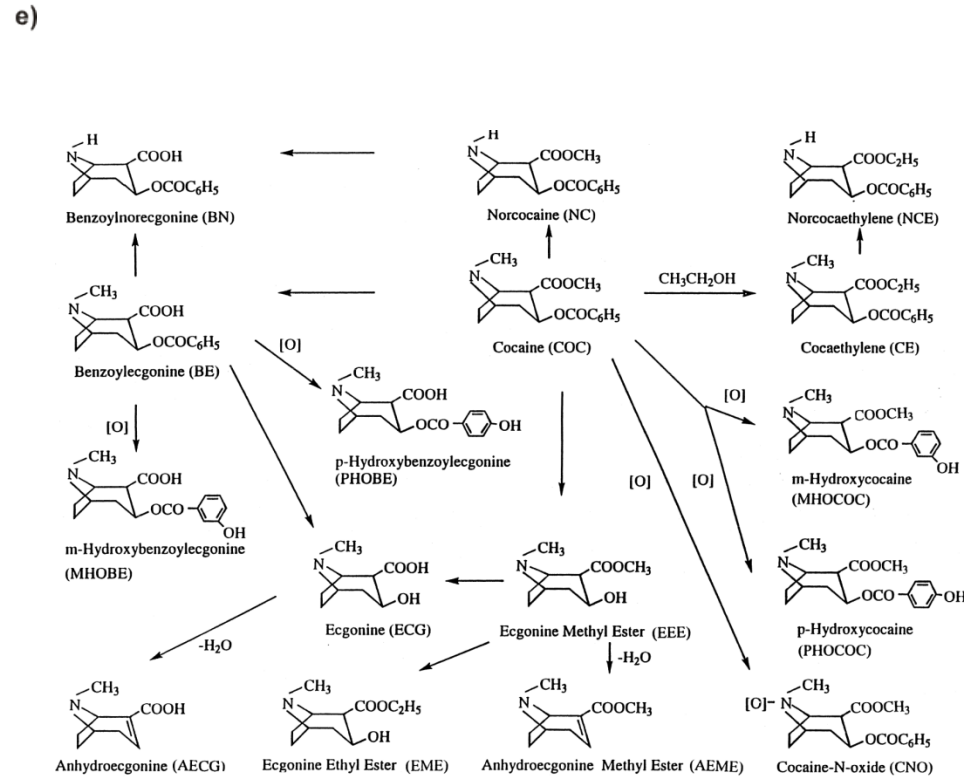
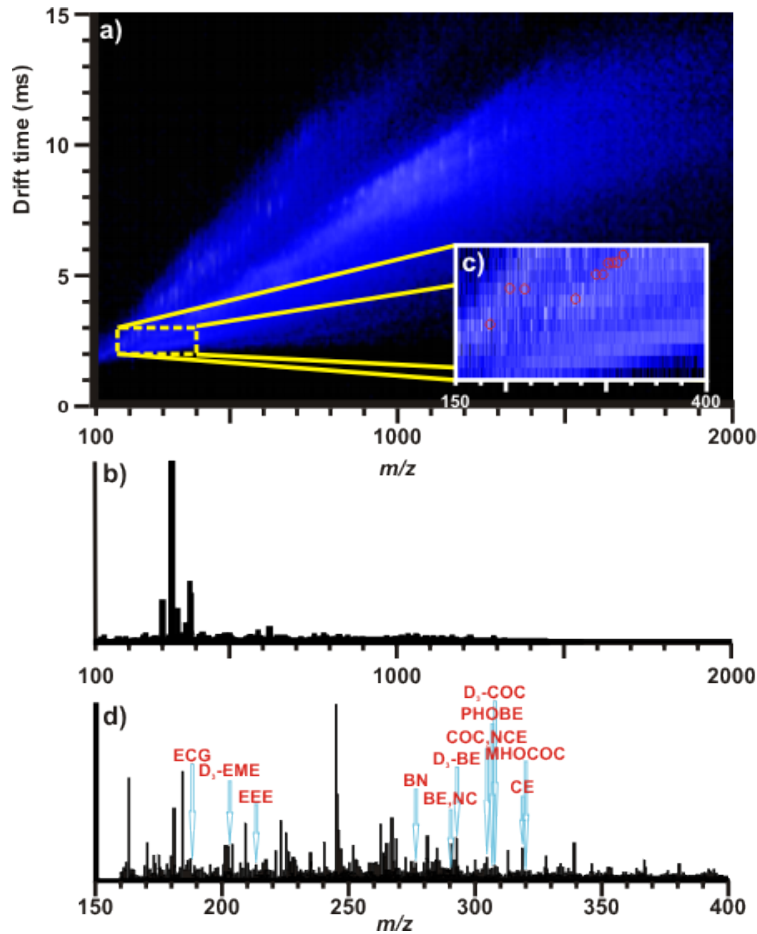
Can leukocytes encode long-term metabolic changes to drug exposure? rat cocaine model



Can leukocytes encode long-term metabolic changes to drug exposure? rat cocaine model

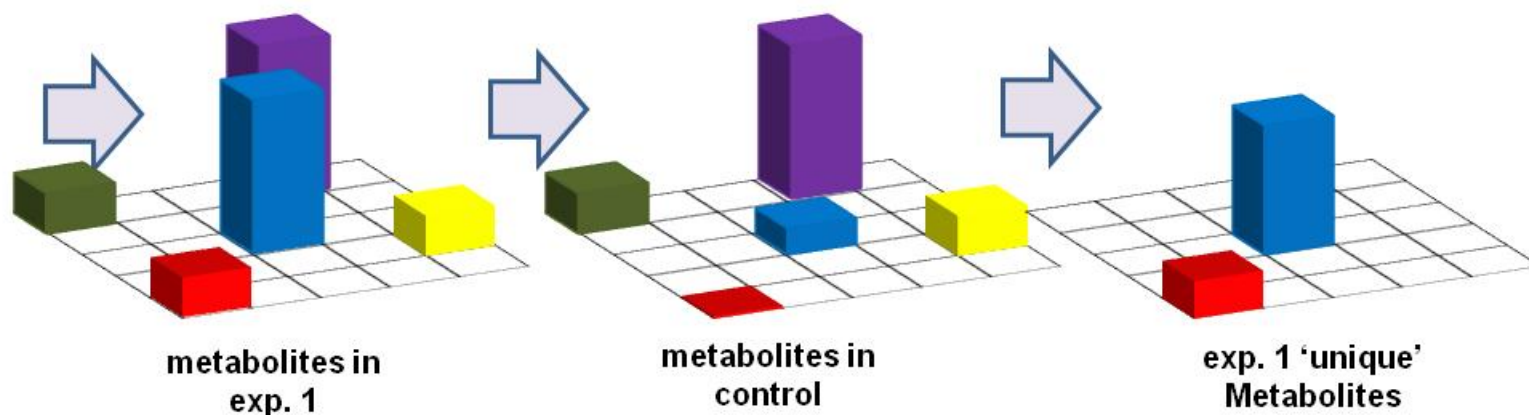
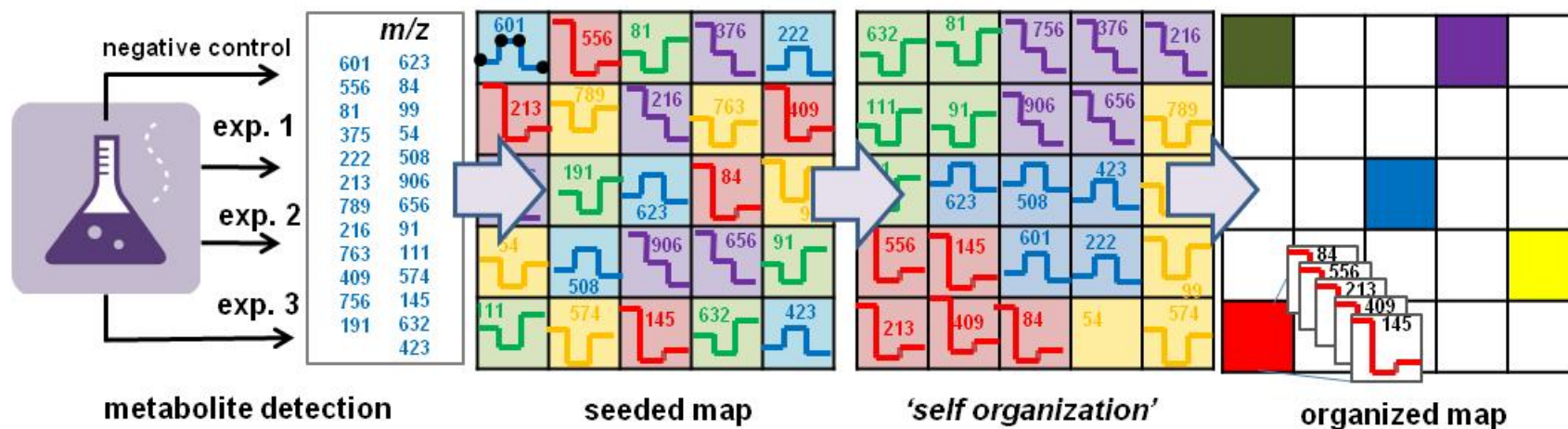


Cocaine metabolism in rat serum

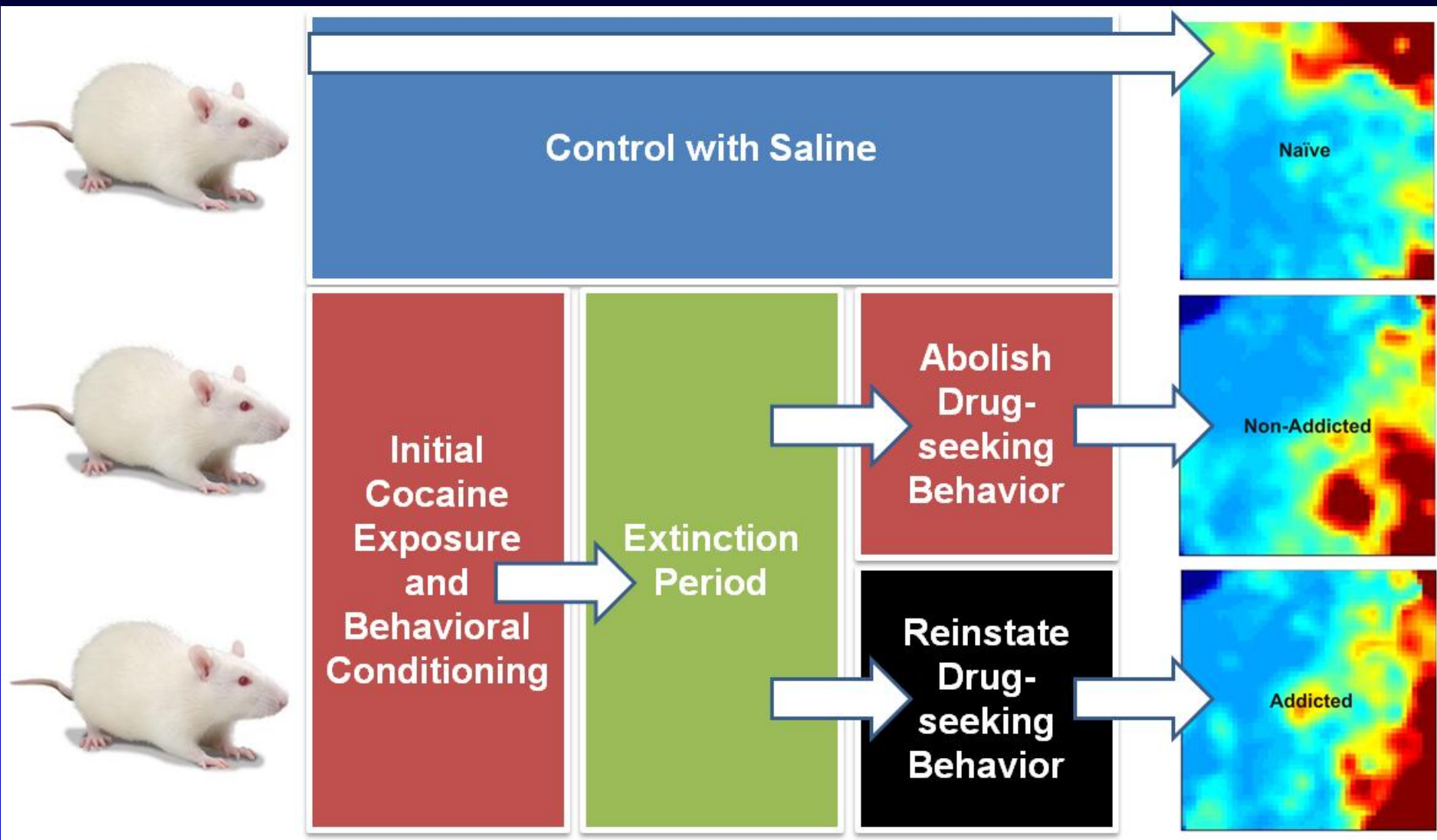


Collaboration with Prof. Shramm-Saypta
(Duke University)

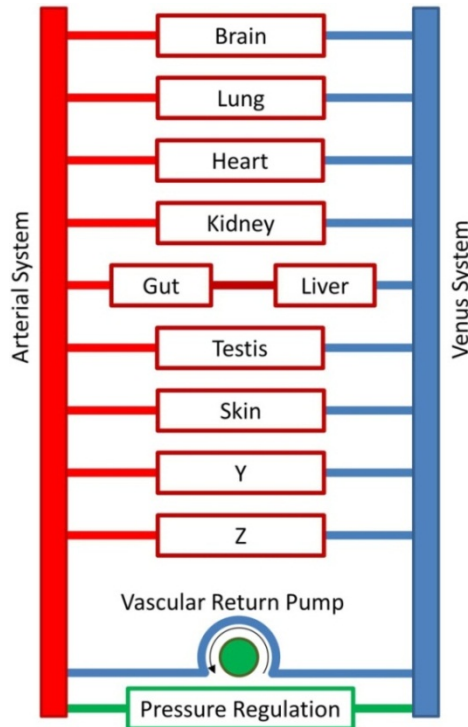
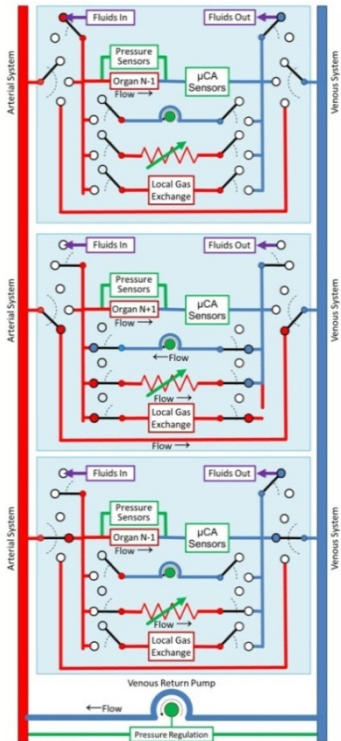
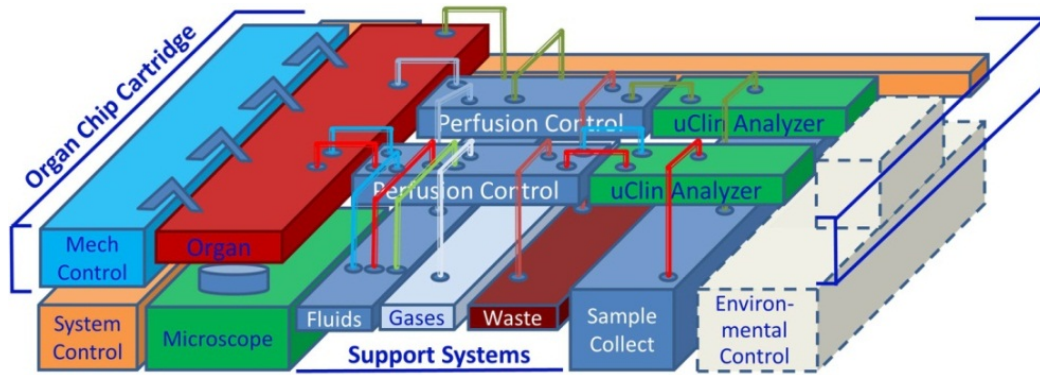
Metabolomic dynamics inspector (MEDI) self-organizing feature maps for target prioritization



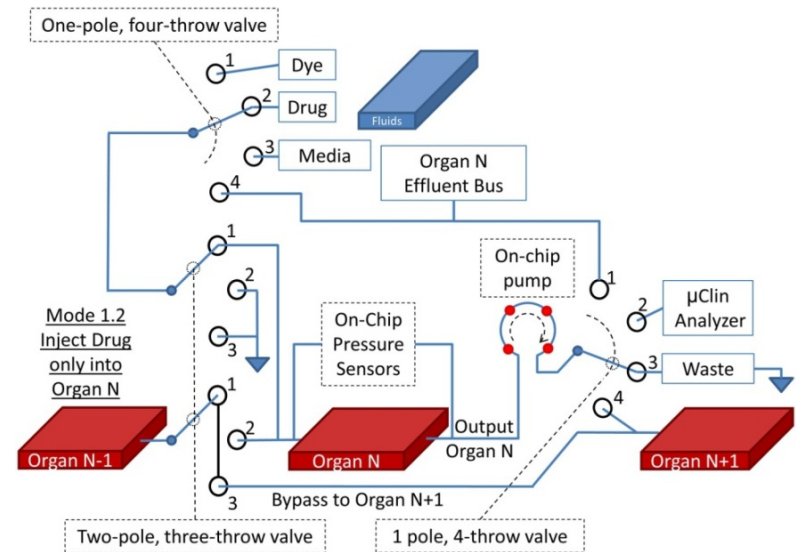
Can leukocytes encode long-term metabolic changes to drug exposure? rat cocaine model



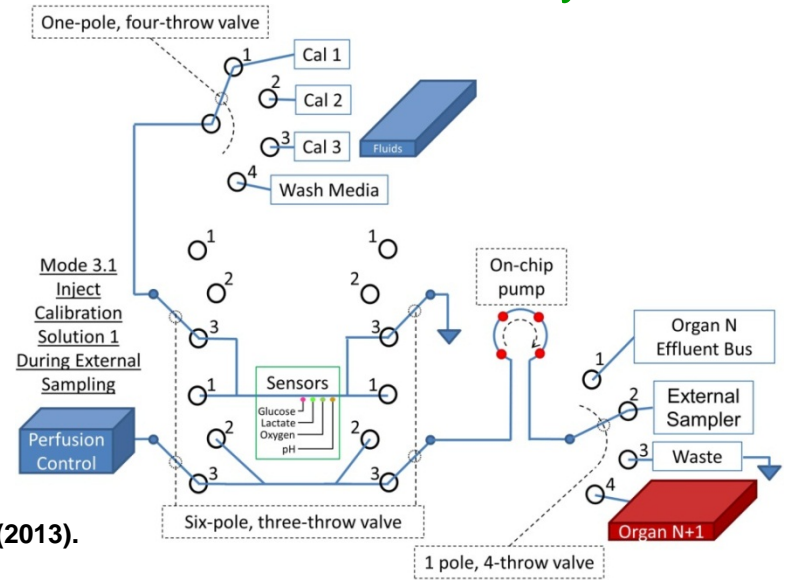
Generalized scheme of the μ Human



Perfusion Controller



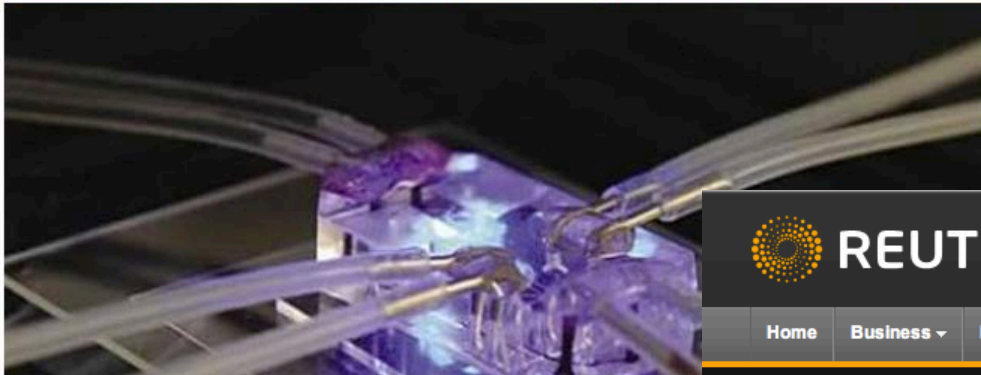
MicroClinical Analyzer



The μ Human and mHuman in the press...

Bloomberg Businessweek Technology

Global Economics Companies & Industries Politics & Policy Technology Markets & Finance Innovation Lifestyle



msn Hotmail More bing TODAY Rock Center Nightly News Meet the Press Dateline Morning J

Home US World Politics Business Sports Entertainment Health

Innovation on NBCNEWS.com

Human Body on a Microchip May Test Drugs



REUTERS

EDITION: U.S.

Register Sign In Search

Home Business Markets World Politics Tech Opinion Breakingviews Money Life Pictures

VIDEO: MOST POPULAR

August 16, 2012



FIND A STATION

home news arts & life music prog

News > Science > Research News

Twitter (9) Facebook (65) Share Comments (2) Recommend (4)

Building Organs, On One Microchip At A Time

DiscoveryNews

EARTH SPACE TECH & GADGETS

Discovery News > Tech News > Scientists Build Living Human Gut on a Chip

SCIENTISTS BUILD LIVING HUMAN GUT ON A CHIP

Analysis by Jesse Emspak
Wed Mar 28, 2012 10:01 AM ET

Artificial jellyfish swims in a heartbeat

Date: Jul 22, 2012

Creation is an amalgam of silicone polymer and heart muscle cells

Chip-based... to revolution... development

June 10 - Researchers are developing microchips of functionality of human organs represent an advance the pharmaceutical companies dollars testing new drugs ineffective - animal trials Transcript)

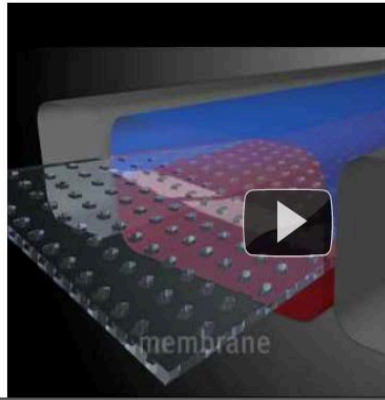
Like 360

MORE VIDEOS

BUSINESS | 7/31/2012 @ 12:50PM | 2,206 views

Military's 'Body-on-a-Chip' Could Fast-Track Pharmaceuticals

+ Comment now

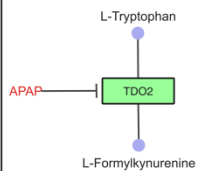
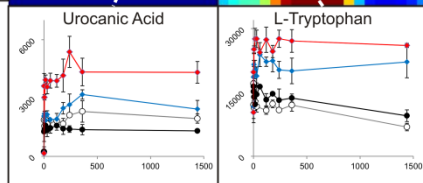
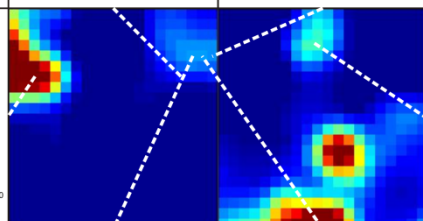
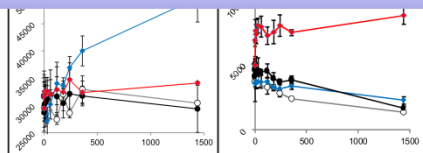


membrane

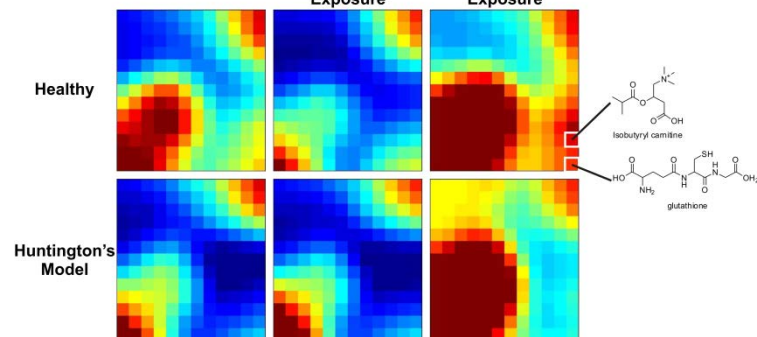
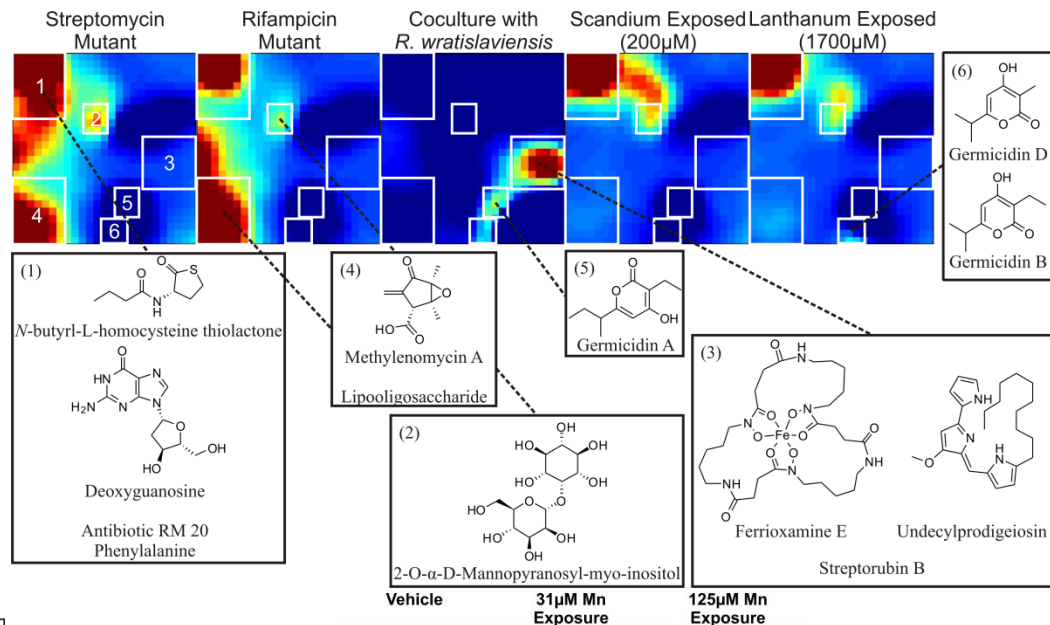
Integrating biology: Translating data to information (Self-organizing maps for untargeted analyses to target identification)

- ❖ Self-organizing maps to elucidate salient molecular features
- ❖ Heat map features direct target species for identification
- ❖ Initially untargeted analyses for panomic molecular coverage

Synthetic biology: liver toxicity to acetaminophen (APAP) by IM-MS



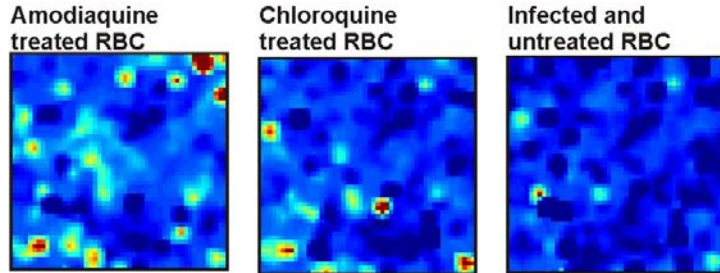
Drug discovery using IM-MS



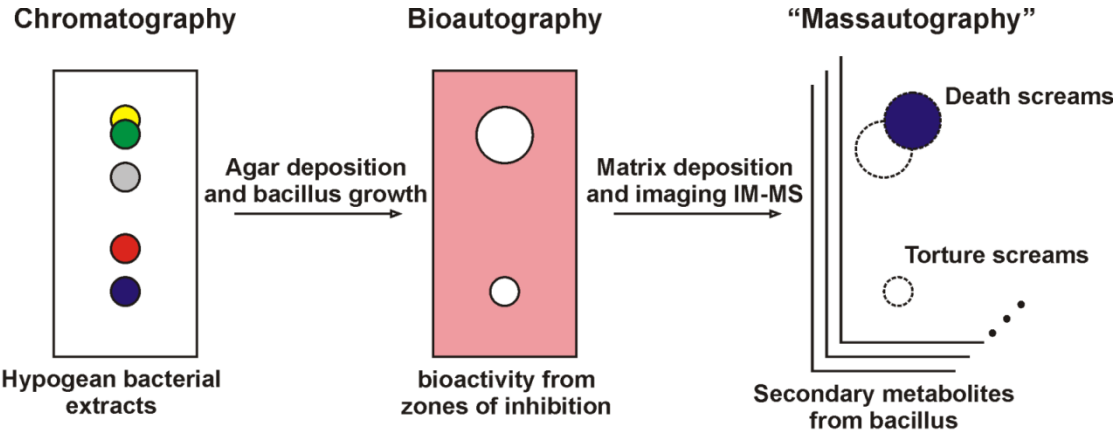
Disease diagnostics with IM-MS

Can we get to systems, synthetic, and chemical biology with integrated omics and informatics?

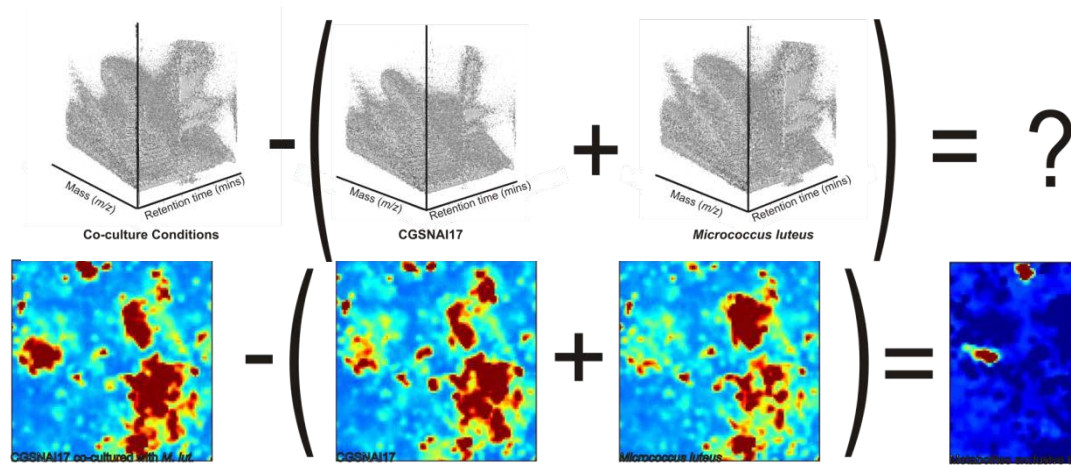
Integrating biology: feature identification



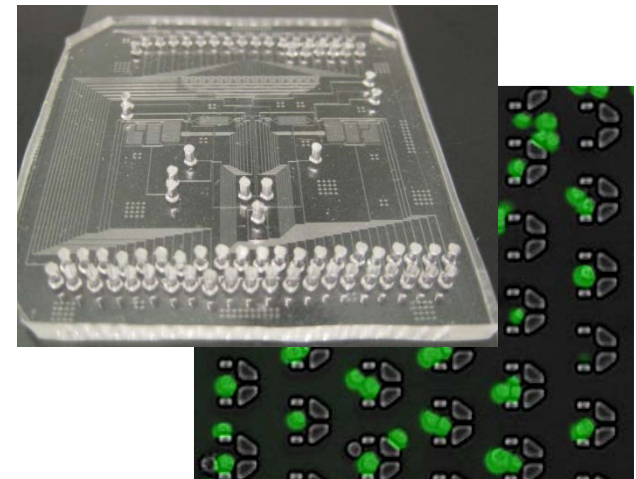
Massautography and imaging mass spectrometry



Drug discovery from complex biological



Metabolic dynamics of cellular perturbation



Laboratory for Structural Mass Spectrometry

Prof. Jody C. May
Dr. Cody R. Goodwin
Ms. Kelly Hines
Mr. Seth Byers
Mr. Jay Forsythe
Ms. Nichole Lareau
Ms. Katie Leaptrot
Mr. Raphael Montenegro
Ms. Sarah Stow
Prof. Sevu Sundarapandian - UHK
Dr. Jeffrey Enders - Ameritox
Dr. Larissa S. Fenn - Waters Inc.
Dr. Randi L. Gant - Aegis Inc.
Dr. Kellen Harkness - EPFL
Dr. Thomas J. Kerr - Optima Inc.
Dr. Michal Kliman - Allergan

Natural product discovery-
Prof. Brian Bachmann, VU Chemistry

Instrumentation-
Dr. John Fjeldsted, Agilent Technologies
Dr. George Stafford, Agilent Technologies
Dr. Ken Imatani, Agilent Technologies
Dr. Ruwan Kurulugama, Agilent Technologies

Computational structural biology-
Prof. Terry Lybrand, VU Chemistry

Dynamic systems biology-
Prof. John Wikswow, VU Physics
Prof. Hod Lipson, Cornell
Dr. Rashi Iyer, LANL
Prof. Don Ingber, Harvard

