

**CLOSE
TO THE
HEART**



FROM VIAL TO ASSAY



Close to the
Heart

Industry Workshop:

Advances in Cardiac Pacing

26th of April 2017

Pluriomics Headquarters

Leiden, The Netherlands

contact: elena@nanion.de

Talk to the experts:



Jim Filippini

sales director, Axion Biosystems, Inc.



Tom O'Brien

CEO, Axion Biosystems, Inc.



Elena Dragicevic

Senior Scientist, Nanion Technologies



Stefan Braam

CEO, Pluriomics

 **pluriomics**

nan]i[on

 **AXION**
BIOSYSTEMS

Lab Demonstrations:

- Maestro,
768- multi-channel MEA
- Pluricyte[®] Cardiomyocytes
- **NEW**
Maestro stimulation plate:
E-Stim MEA 48-well plate

Assessment of pro-arrhythmic effects using
Pluricyte[®] Cardiomyocytes on the
Axion Maestro system

