

nanjion

# What's New in CiPA: An Update of Stem Cell and Ion Channel Assays for Safety Studies



## Nanion Hosted Session:

**Bernard Fermini, Pfizer:**  
"The CiPA Ion Channel Working Group: Current status and next steps."

**James Kramer, Charles River Laboratories:**  
"Challenges for Pharmaceutical Industry and Opportunities for Safety Pharmacology"

**Herbert Himmel, Bayer Pharma:**  
"From ICHS7B to CiPA – the changing landscape of cardiac ion channel assessment"

**Chairs: Sonja Stölzle-Feix, Corina Bot, Atsushi Ohtsuki, Nanion Technologies**  
**Monday 19<sup>th</sup> of September,**  
**18:00 – 19:00 pm, room 7**

**Sonja Stölzle-Feix, Nanion":**  
"Cardiotoxicity screening with Cellartis cardiomyocytes: combining electrophysiology with impedance"

**Monday 19<sup>th</sup> of September,**  
**7:00–8:00 am, room 16**

## Posters:

**# 0129: An "All Inclusive" Package for Cardiac Safety: The Six Big on One Automated Patch Clamp Chip**

**# 0128: Impedance and Extracellular Field Potential for Cardiac Safety Assays: A Combined Approach for NonInvasive Screening of iPS Cells**

**# 0242: Human ventricular stem cell cardiomyocytes: In vitro assays and screening platforms for pro-arrhythmia risk prediction**



Don't miss:

Nanion's Hosted Company Session  
Nanion's Poster on cardiac safety  
Nanion's Networking Reception

**Takara Clontech  
Hosted Session:**