



nanjion

STEM CELL THEMED CONFERENCE

8th of March 2017

Granta Park, Cambridgeshire

Contact: elena.dragicevic@nanion.de

Presenter:

- Mohamed Kreir (Janssen)
- Tristan Prichard-Meaker (Axiogenesis)
- Michela Nosedà (Imperial College London)
- Sonja Stölzle-Feix (Nanion Technologies)
- Matthew Bridgland Taylor (AstraZeneca)
- Said El Haou (Metrion Biosciences)

Lab Demonstrations:

- Patchliner
- CardioExcyte 96
- Maestro