

# CardioExcyte 96 SOL

Optogenetics meets cardiac safety



This is  
New:

Optical pacing

Add-On for the  
CardioExcyte 96

long-term  
stimulation during  
maturation cycles

Visit us at booth #405  
to see our new  
CardioExcyte 96 SOL

**The new CardioExcyte 96 SOL** uses LEDs for spatially uniform stimulation of cells transfected with light-gated ion channels such as Channelrhodopsin2 (ChR2).

**CardioExcyte 96 SOL** offers you:

- Mechanistic insights into contractility and electrophysiology of iPS cardiomyocytes
- Pacing without stimulus artifacts
- Pacing without electrical stress for cells

nanjion

**Nanon Europe**  
info@nanion.de  
Phone: +49 89 2 18 99 79 72  
www.nanion.de

**Nanon USA**  
info@naniontech.com  
Phone: 1-888-9-NANION  
www.naniontech.com

**Nanon China**  
info@nanion.cn  
Phone: +86 10 82 17 63 86  
www.nanion.cn

**nanjion**

# Optical stimulation and investigation of impedance and extracellular field potentials in parallel

*"The CardioExcyte 96 is a hybrid system capable of recording impedance and extracellular field potentials of beating networks of cells. The advantages of optical stimulation over electrical stimulation include the highly precise timing, all cells are stimulated exactly at the time of the trigger. In contrast, electrical stimulation propagates from the electrode across the well and thus cells are stimulated progressively. A mean beat*

*calculation of precisely timed beats enables in - depth compound analysis and concentration response dependencies which is one of the key software capabilities of the CardioExcyte Control/ DataControl software package that is provided with each CardioExcyte 96 system."*

**Dr. Sonja Stoelzle Feix, Director Scientific Affairs, Nanion Technologies GmbH**

## Do not miss:

- To learn about the advantages of optical stimulation in comparison to electrical stimulation of cardiomyocytes
- To learn about frequency-dependent compound effects

## Visit our Exhibitor presentation (Monday, 12:30 PM, room 221):

- "Ion channel Drug Discovery: Beyond the bottleneck and ready for CiPA"

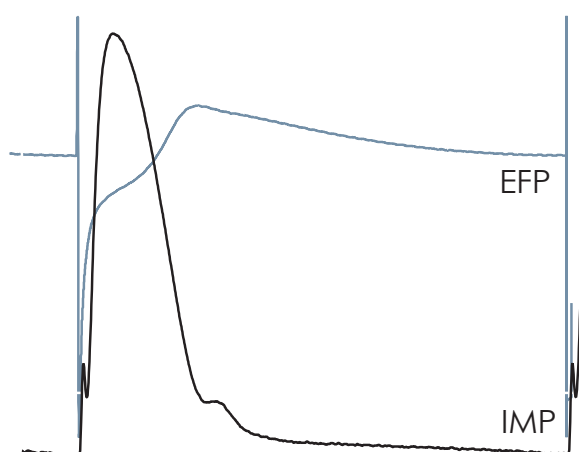
## Visit our booth # 405 and talk to our CardioExcyte 96 specialists:

- Corina Bot, Niels Fertig, Michael George

## Visit our platform session oral presentation (Monday, 9:45 AM, room R04/05)

- Corina Bot: "Optical stimulation of iPS cardiomyocytes allows brand new insights into contractility and electrophysiology conjunctions" (Program # 797-Plat)

Timed stimuli allow exact IMP and EFP mean beat overlays of the same cells



Nifedipine (black) shorten FPD and reduce contraction strength, 2 Hz paced cells

