

# Label-free analysis of Na<sup>+</sup>/Ca<sup>2+</sup>-exchanger (NCX) isolated from iPSC-derived cardiomyocytes



Maria Barthmes, Andre Bazzone, Ulrich Thomas, Andrea Brüggemann, Michael George, Niels Fertig, [Alison Obergrussberger](#)

Nanion Technologies GmbH, Ganghoferstr. 70A, 80339 Munich, Germany, contact: [Info@nanion.de](mailto:Info@nanion.de)

## Abstract

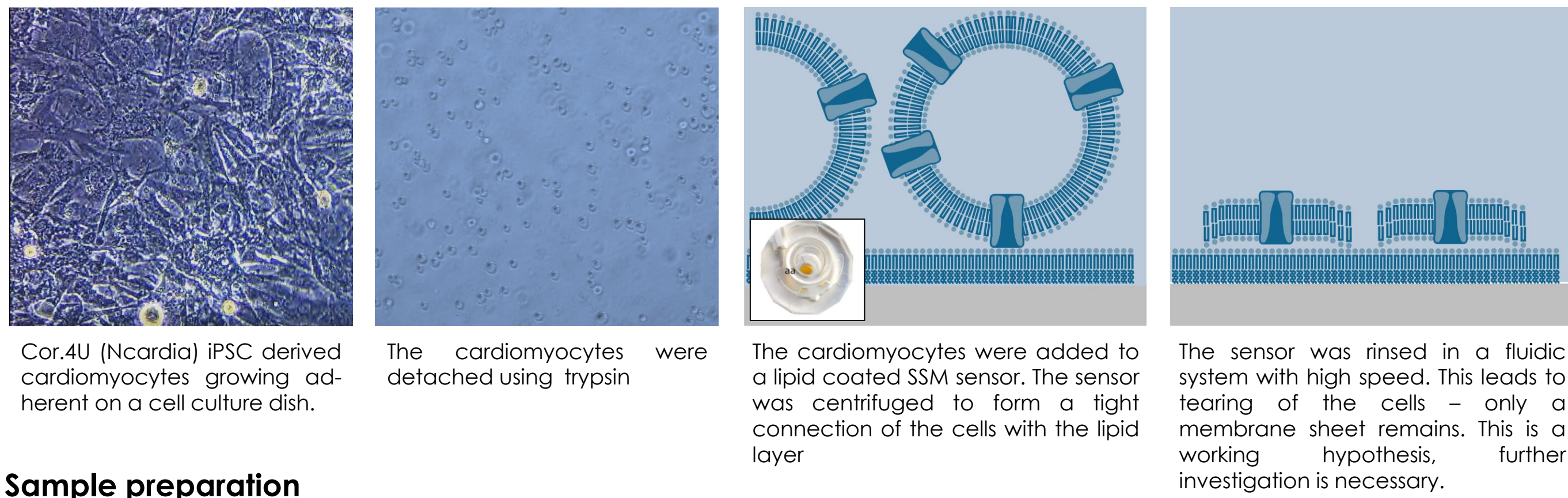
The Sodium-Calcium Exchangers (NCX) play an important role in the cellular calcium homeostasis under physiological and pathological conditions. NCX has been of interest as a pharmacological target for many years, in particular because clinical trials involving inhibitors of the sodium-proton exchanger, NHE, have delivered mixed results. Inhibition of the reversed mode of NCX is thought to be beneficial in ischemia/reperfusion injury by reducing cardiac, neuronal and renal infarct areas. Moreover, inhibition of NCX has been proposed to exhibit an anti-arrhythmic effect and therefore, may provide a novel target for the treatment of a variety of arrhythmic pathologies. So far, a number of studies have shown promising results but investigations are limited by the currently available NCX inhibitors such as KB-R7943, SEA-0400 and SN-6 which are only partially specific.

To drive the progress in pharmacological NCX research, new methods to measure NCX function are needed. At the current time, functional investigation of NCX range from patch-clamp, calcium flux assays, Langendorff-perfused hearts to studies in whole animals. We have developed an electrophysiological method to investigate NCX function which is based on the solid supported membrane (SSM) technology. HEK cells overexpressing NCX or human induced pluripotent stem cell-derived cardiomyocytes (hiPSC-CMs) were used on a single-well or 96-well SSM electrophysiology device and NCX was recorded from these cells. NCX was activated using Ca<sup>2+</sup> in the buffer and inhibited by Cd<sup>2+</sup> and other compounds.

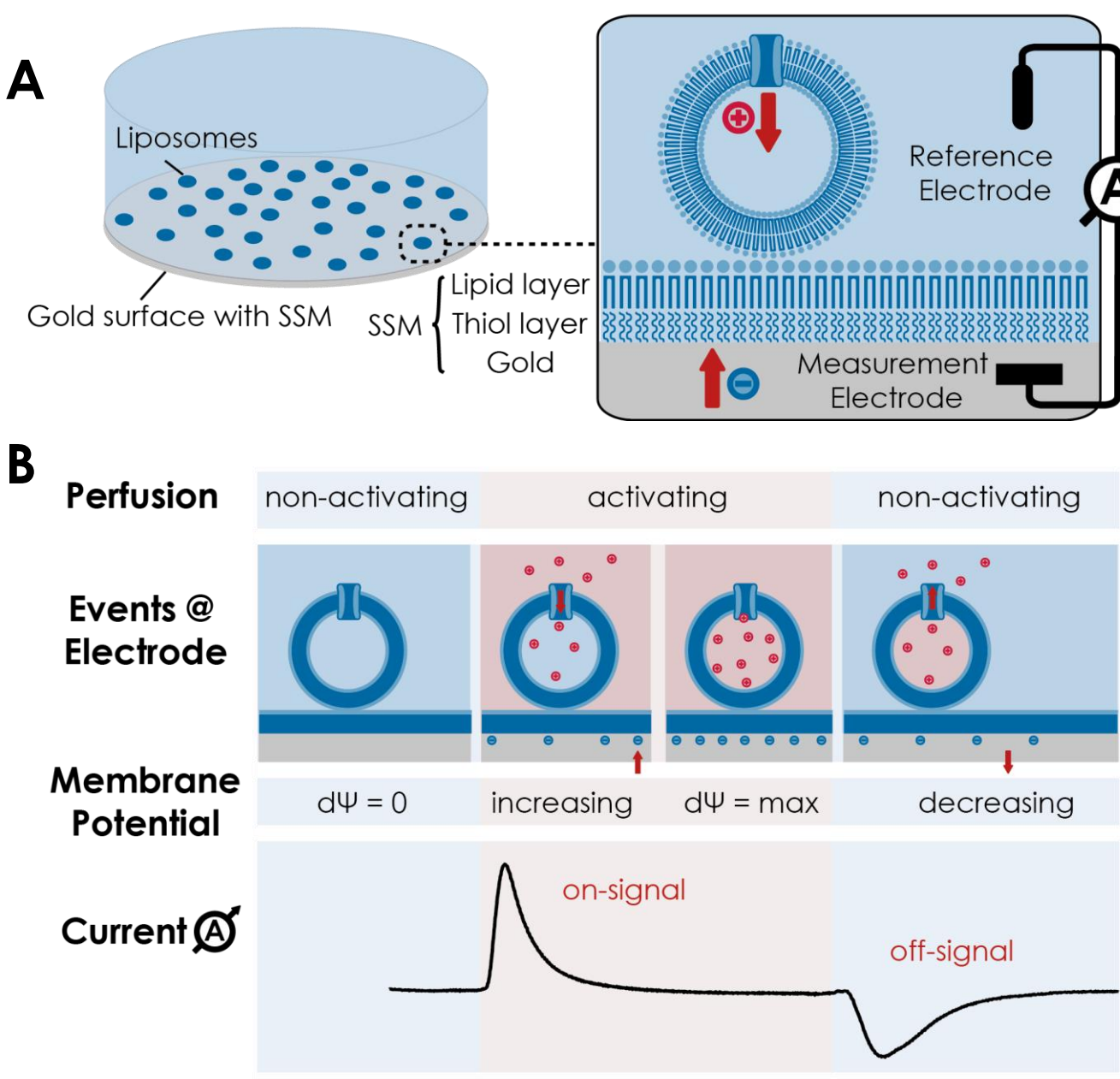
## Measuring NCX activity in human iPSC derived cardiomyocytes

Human iPSC derived cardiomyocytes are being investigated as a model for cardiac safety assessment. To measure native NCX in these cardiomyocytes a cell based assay was developed. Cardiomyocytes were detached from the culture dish and added to the lipid coated SSM sensor. Where the cell connected with the lipid layer. Upon rinsing of the sensor

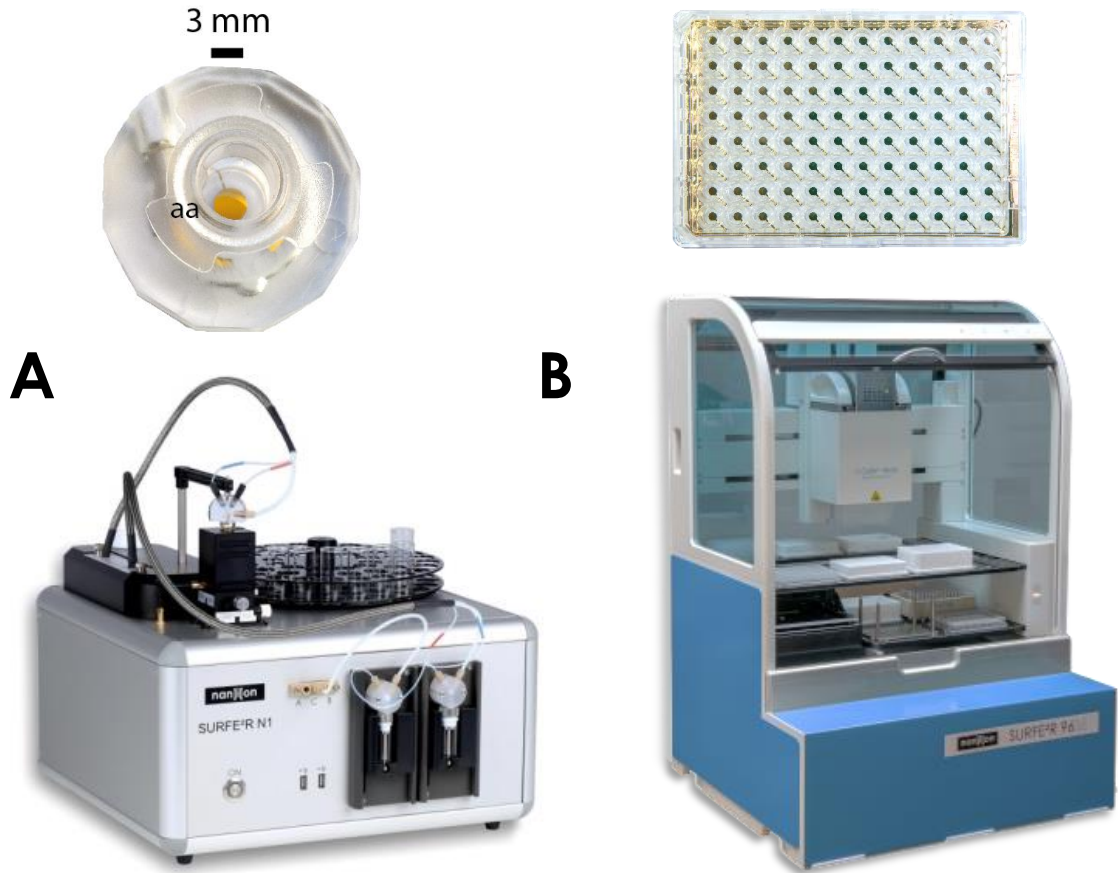
with high fluidic speed, the cells detached again, but a sheet of the cell membrane remains on the sensor. NCX currents can be evoked in these sheets. For a higher NCX signal female cardiomyocytes were used. This method enables the efficient investigation of the isolated cardiac NCX current in a native membrane.



## Resolving low amplitude ion currents



**"Solid Supported Membrane (SSM)-based electrophysiology"** is a method, which allows the resolution of low amplitude electric events in biological membranes, bringing the advantages of electrophysiology to the transporter field. It enables real-time activity measurement of electrogenic SLC-transporters, MFS-transporters and ion pumps, localized in the plasma membrane, in intracellular or bacterial membranes. Unlike cell-based electrophysiology, a large capacitive sensor coated with a high amount of membrane vesicles or proteoliposomes is used. In this way any process that generates an electrical potential can be registered with high amplification. The method was established in the 90's and has only recently been scaled-up to a 96-well format.



Instrumentation: We used the SURFER N1 (A) and the SURFER 96SE (B) instruments in this project. The respective SSM sensors are depicted above the instruments. The SURFER N1 is a flexible single channel measurement instrument, whereas the SURFER 96SE is a robotic device designed for experiments that require higher throughput. C: Membrane proteins successfully studied using SSM-based electrophysiology

Transporters			
Inorganic Ions	Amino acids	Sugars	Organic Ions
NhaA	PEPT1 (Slc15a1)	SGLT1/2 (Slc5a1/a2)	OC12 (Slc22a2)
NhaP	YdR	MeIB	CNT1 (Slc28a1)
NhaB	YhP/DtpB	LacY	ANT (Slc25a4)
NCX1 (Slc8a1)	PuP	FucP	AAC
Cic-7	GHF	XyIE	GAT1 (Slc6a1)
EcCic	EAAC1 (Slc1a1)	GlcP	BeP
NrC	PAT1 (Slc36a1)		CHT (Slc5a7)
Am11-3	ArcD		NupC
Am1B	CAZB (Slc7a2)		NacT (Slc13a5)
SuIP	GlyT1/2 (Slc6a7/a5)		
NIS (Slc5a5)			
NaP2b (Slc34a2)			
MnH2			

ATPases		Redox-driven ion pumps	
NaK-ATPase		Complex I	
HK-ATPase		respiratory chain complex I/II	
SERCA		respiratory chain complex II/III	
V-ATPase		cytochrome c-oxidase	
FoF1-ATPase		respiratory chain complexes I/II/IV	
Kdp-ATPase			
CopA			
ATP7A/B			
VrPPase			

Channels and Pores	
Gramicidine	
P2X2	
nACHR	
A/M2	
UCP1 (Slc25a7)	
TRPC5	
TRPA1	
CFTR	
AQP6	

Light-driven ion pumps	
Bacteriorhodopsin (BR)	
Oxymis marina Rhodopsin	
Rhodopsin-2 (KR2)	
Halorhodopsin (HR)	
Acerhodopsin	
Channelrhodopsin (ChR)	

## Summary and Future prospects

We have developed two novel methodic approaches for the functional investigation of NCX and have recorded NCX in HEK cells and hiPSCs. NCX was activated by Ca<sup>2+</sup> in the buffer and could be inhibited by a number of pharmacological agents.

### Future Experiments

- Further validation and optimization of the methods is required.
- We intend to develop a third complementary method using a MEA-like system to investigate the influence of NCX inhibitors and cross reactions of NCX inhibitors and pro-arrhythmic drugs on the beating behavior of iPSC derived cardiomyocytes.

- The developed methods shall be applied in a project to characterize drug effects on NCX in more detail.
- The next step intended is to screen the proarrhythmic drugs of the CiPA initiative compound list using the high throughput method for effects on NCX.

### Contact us if you are interested in collaborating with us to:

- Apply the high throughput method for the development of novel NCX inhibitors.
- Evaluate the need for NCX safety testing of new drug candidates using the cardiomyocyte based method.

## Literature

Amran et al., Cardiovascular Drug Reviews, Vol. 21, No. 4, pp. 255-276  
Antoons et al., Pharmacology & Therapeutics 134 (2012) 26-42  
Hermann et al., Cardiovascular Research (2013) 99, 780-788  
Huang, JPTM 87 (2017) 11-23  
Lee and Hryshko, Cardiovascular Drug Reviews, Vol. 22, No. 4, pp. 334-347  
Pignataro, Translational Stroke Research, DOI 10.1007/s12975-013-0308-8  
Watanabe, et al., J. Pharmacol.Sci.102, 7 - 16 (2006)  
Matsuda, et al., 2001, JPET, 298(1): 249-256  
Watanabe & Kimura., 2001, Jpn. J. Pharmacol. 85: 370-375

## Development of a screening tool for NCX

To screen compounds for an effect on NCX for safety purposes or during development of novel inhibitors a robust method with a high throughput is required. We employed the 96-well based platform SURFER 96SE to record NCX activity from stably transfected HEK293 cells. The cells were cultured, harvested and a membrane preparation was performed. This has the advantage that the resulting signal amplitudes are higher, no running cell culture is required and it yields a very high reproducibility. The membrane preparation is defrosted and added to a 96-well sensor where it is recorded by fluidic activation with calcium.

