



Symposium & Wet Lab Demonstrations

Nanion 7th Annual User Meeting

Venue:

- ⇒ The 7th Annual User Meeting will be held at Nanion Headquarters, Gabrielenstrasse 9, 80636 Munich, Germany
- ⇒ October 2016: Wednesday the 26th (1:00 pm) – Thursday the 27th (2:00 pm)

Preliminary program:

Symposium:

- ⇒ **Gerard Baaken**, Ionera: "Orbit16/Mini: Generic tools for the analysis of pore forming proteins"
- ⇒ **Matt Burnham**, AstraZeneca: "Towards exploration of 3D iPSC-CM models with the CardioExcyte96."
- ⇒ **Teun de Boer**, Univ. Utrecht: "Dynamic clamping on an APC machine: adding virtual IK1 channels to cardiomyocytes."
- ⇒ **Divya Mirrington**, Univ. Nottingham: "Modelling the cardiac channelopathies using human induced pluripotent stem cells"
- ⇒ **Fabian Haeusermann**, Roche (topic: CardioExcyte)
- ⇒ **Chris Mathes**, Icagen: "Characterization of subtype selective KCNQ activators using the Syncropatch 384PE."
- ⇒ **Stephen Tucker**, Univ. Oxford: "Port-a-Patch Recordings of Purified TREK-2 K⁺ Channels Reveal an Asymmetric Mechanosensitivity"

Wet Lab Demonstrations:

- ⇒ Port-a-Patch, Patchliner, CardioExcyte 96, SyncroPatch 384PE, Orbit Mini, SURF2R SE, Maestro



Symposium & Wet Lab Demonstrations

Accommodations:

- ⇒ Hotel Orly (Gabrielenstrasse 6, 80636 Munich)
<http://www.hotel-orly.de>
2 minutes walking distance from Nanion headquarters
Single room: 105,00 €, double room: 110,00 € (breakfast included)
- ⇒ Novotel München City Arnulfpark (Arnulfstrasse 57, 80636 Munich)
<http://www.novotel.com/de/hotel-8866-novotel-muenchen-city-arnulfpark/index.shtml>
10 minutes walking distance from Nanion headquarters
Single room: 124,00 €, double room: 143,00 € (breakfast included)

Travel:

- ⇒ If you travel by **train** (ICE): Either change to the „S-Bahn“ at Munich main station and exit at „Donnersberger Brücke“ (2 stations) or change to the „S-Bahn“ in Pasing and exit at the „Donnersberger Brücke“ (3 stations). From here, you can reach both hotels and nanion headquarters within a walking distance of max.10 minutes.
- ⇒ If you travel by **airplane**: Change at the airport Munich to S-Bahn „S8“ and exit at „Donnersberger Brücke“ (approx. 75 minutes)

Poster:

- ⇒ If you would like to present a poster (DINA0 portrait orientation), please send a title and an abstract to: gesa.rascher-eggstein@nanion.de

Oral Presentation:

- ⇒ If you would like to give an oral presentation, please forward a title and an abstract to: gesa.rascher-eggstein@nanion.de



Symposium & Wet Lab Demonstrations

Co-marketing opportunities for collaboration partners:

- ⇒ If you like to bring one **rollup banner**, it will be placed within the main room.
- ⇒ If you like to present a **poster**: We are delighted to accept poster, preferably with data on nanion technology or helpful content for our attendees.
- ⇒ If you like to bring **marketing material** as flyer and brochures: We will prepare one table for all collaboration partners on which marketing material can be placed.

If you are interested in any of the above co-marketing opportunities, please inform me in advance via e-mail to ensure an optimal organization of the User Meeting: gesa.rascher-eggstein@nanion.de

- ⇒ Collaboration partners: Please send us a **company description** with max. 100 words and your **logo**, as we would like to implement both in our User Meeting booklet.

If you have any questions, I am happy to help.

Please do not hesitate to call me directly or send me an e-mail:

- ⇒ phone: +49 89 2189979-73
- ⇒ mobile: +49 176 396-08117
- ⇒ e-mail: gesa.rascher-eggstein@nanion.de

Looking very much forward to seeing you end of October!

Gesa Rascher-Eggstein

Marketing Manager Nanion Technologies GmbH