

Satellite Symposium Announcement

New and Emerging Techniques in Electrophysiology

Sponsored by: Nanion Technologies GmbH, Munich
and the EU-project MRTN-CT 2005-019335 Translocation

Saturday 25.03.2006, 10:00-evening

Physiology Dept.,
Pettenkoferstr. 12,
University Munich

Co-Chairs: Niels Fertig (Munich), Jan Behrends (Freiburg)

Program:

10:00-10:05
Welcome Address

10:05-11:00
Richard Wagner (Osnabrück), Keynote

"Simultaneous optical and electrical recordings from nano- μ m scaled horizontal bilayers."

11:00-11:30
Rikard Blunck (UCLA)
„Fluorescence measurements from planar lipid bilayers“

11:30-12:00
Markus Sondermann (Freiburg)
„Patch clamp on a chip: single channel experiments
in solvent-free lipid bilayers“

12:00-13:00
Lunch Break

13:00-13:30
Stefan Huber (Tübingen)
„Electrophysiology of malaria-infected erythrocytes“

13:30-14:00
Gregor Dernick (Basel)
„Patch Amperometry“

14:00-14:30
Jörg Geiger (Frankfurt)
„Patch clamp recordings from a synaptic bouton“

14:30-15:00
Coffee Break

15:00-15:30

Clemens Möller (Hamburg)

„Ion Channel Assays in Drug Discovery and Development”

15:30-16:00

Michael Fejtl (Tübingen)

„ChipTips allow fast solution exchange for the Flip-the-Tip technology”

16:00-16:30

Bela Kelely (Frankfurt)

„SURFE²R technology - Cell free electrophysiology for the study of transporters”

there will be a small lunch snack and in the evening a
Brez'n & Beer reception
(featuring a great Bavarian Monastery Brew (dark))

if you plan to attend, please briefly send an email with subject “symposium” to **info@nanion.de**