

Nanion-TWMU Ion Channel Forum 2017

- Automated Patch Clamp and beyond -

Date : Thursday, 27.7.2017, 10:00~17:00

- Reception dinner will take place at the KEIO PLAZA HOTEL TOKYO, Shinjyuku
- E-Physis Workshop will be held on Friday, 2017.7.28

Location : ABMES Conference Room(02N501), TWIns 2F
Institute of Advanced Biomedical Engineering and Science
Tokyo Women's Medical University
2-2 Wakamatsu-cho, Shinjyuku-ku, Tokyo 162-0056

Capacity : 100 seats



Confirmed Speakers:

Dr. Ryo Matsumoto

Pharmacological Group, Pharmaceutical Research Department, Biological Research Laboratories, NISSAN CHEMICAL INDUSTRIES. LTD.

Katsuhisa Matsuura, M.D., Ph.D.

Institute of advanced Biomedical engineering and sciences, Tokyo Woman's Medical University

Sayo Suzuki, Ph.D.

Division of Pharmaceutical Care Sciences, Center for Social Pharmacy and Pharmaceutical Care Sciences,

Keio University

Dr. Ayumi Hirano-Iwata

Advanced Institute for Materials Research(AIMR), Tohoku University

Reserch Institute of Electrical Communication, Tohoku University

Toshikazu Nishie, Ph.D.

CMD Center, TAKARA BIO INC.

Dr. Takuji Yamaguchi

Charles River Laboratories Japan, Inc.

Dr. Niels Fertig,

Founder and CEO Nanion Technologies, Nanion Technologies GmbH

Dr. Sonja Stoelzle-Feix,

Director, Scientific Affairs, Nanion Technologies GmbH

And more !

Email to ohsuki@nanion.jp for your registration !

nan]i[on