

# SOPHION BIOSCIENCE SEMINAR

11  
October  
2019

Venue: Knowledge Capital Conference Rooms, Room C04  
Grand Front Osaka, North Bld., Tower C, 8F  
3-1 Ofuka-cho, Kita-ku, Osaka 530-0011, Japan

## Agenda

10:00-10:30	Registration
10:30-10:40	Welcome remarks and safety briefing Dr. Yuji Tsurubuchi, Country Manager, Sophion Bioscience K.K., Japan
10:40-11:00	TBD Mr. Thais Johansen, CEO, Sophion Bioscience A/S, Denmark
11:00-12:00	Educational lecture: TBD Prof. Fumihito Ono, Department of Physiology, Division of Life Sciences, Faculty of Medicine, Osaka Medical College, Japan
12:00-13:00	Lunch break (Lunch boxes served)
13:00-13:40	Ten years of automated patch-clamp platforms use at Sanofi (Tentative) Dr. Michel Partiseti, Sanofi, France
13:40-14:20	TBD Dr. Edward Humphries, Metrion Biosciences Ltd., UK
14:20-15:00	Updates on QPatch and Qube applications and their new features (tentative) Application Scientist or Product Manager, Sophion Bioscience A/S, Denmark
15:00-15:30	Coffee break (Beverages and sweet treats served)
15:30-16:30	Basic course of training in QPatch/Qube operations and data analyses for new users Dr. Kazuya Tsurudome, Senior Application Scientist, Sophion Bioscience K.K., Japan
16.30-17.00	Tricks and tips for advanced users to maximize utilization of QPatch and Qube Dr. Hironori Ohshiro, Application Scientist, Sophion Bioscience K.K., Japan
17:00-17:10	Closing remarks and adjourn Mr. Naoto Ueda, President & CEO, Physio-Tech Co., Ltd., Japan
17:30-20:00	Reception party at the next-door Room C03 (Beverages and foods served)