

SOPHION BIOSCIENCE SEMINAR

19
October
2018

Venue: CMIC Pharma Science Co., Ltd., Tokyo Office

Hamamatsucho Bldg., 22nd Floor, 1-1-1 Shibaura, Minato-ku, Tokyo 105-0023, Japan

Agenda

09.30-10.00	Registration
10.00-10.10	Welcome remarks and safety briefing Dr. Yuji Tsurubuchi, Country Manager, Sophion Bioscience K.K., Japan
10.10-10.30	Sophion Mr. Thais Johansen, CEO, Sophion Bioscience A/S, Denmark
10.30-10.50	Corporate Outline (Tentative) Dr. Yasuki Akie, CMIC Pharma Science Co., Ltd., Japan
10.50-12.00	Educational lecture: Stoichiometry and functions of potassium channel auxiliary subunits Prof. Koichi Nakajo, Division of Integrative Physiology, Department of Physiology, Jichi Medical University, Japan
12.00-13.00	Lunch break (lunch boxes served)
13.00-13.40	Utilising the Qube in a CRO environment: Case studies from HTS and hit-to-lead projects Dr. Juha Kammonen, Principal Scientist, Charles River Laboratories, United Kingdom
13.40-14.10	Current and future research & development activities at Sophion Dr. Thomas Binzer, VP, Sophion Bioscience A/S, Denmark
14.10-14.40	Introducing QPatch II / CiPA hERG Milnes kinetic assay on QPatch Dr. Kazuya Tsurudome, Sophion Bioscience K.K., Japan
14.40-15.20	Coffee break (Beverages served)
15.20-16.00	Myocardial contraction assay using imaging technology -SI8000 Cell Motion Imaging System- Dr. Tomohiro Hayakawa, Sony Imaging Products & Solutions Inc., Japan
16.00-16.40	High-resolution <i>in vitro</i> electrophysiology system for detailed characterization of neuronal cells and networks Dr. Marie Engelene J.Obien, VP Marketing and Sales, Maxwell Biosystems AG, Switzerland
16.40-17.00	Closing remarks Mr. Hiroshi Kosaku, President, CMIC Pharma Science Co., Ltd., Japan
18.00-20.00	Reception party (Beverages and foods served)

SOPHION BIOSCIENCE SEMINAR

19
October
2018

会 場: シミックファーマサイエンス株式会社東京オフィス会議室
浜松町ビルディング 22 階 (旧東芝ビル) 東京都港区芝浦 1-1-1

Agenda

09.30-10.00	受付
10.00-10.10	開会の挨拶 ソフィオンバイオサイエンス株式会社 代表取締役 鶴淵 雄士
10.10-10.30	Sophion Mr. Thais Johansen, CEO, Sophion Bioscience A/S, Denmark
10.30-10.50	会社案内 シミックファーマサイエンス株式会社 秋江 靖樹
10.50-12.00	教育講演: Stoichiometry and functions of potassium channel auxiliary subunits 自治医科大学 生理学講座統合生理学部門 中條 浩一 教授
12.00-13.00	昼食
13.00-13.40	Utilising the Qube in a CRO environment: Case studies from HTS and hit-to-lead projects Dr. Juha Kammonen, Principal Scientist, Charles River Laboratories,
13.40-14.10	Current and future research & development activities at Sophion Dr. Thomas Binzer, VP, Sophion Bioscience A/S, Denmark
14.10-14.40	QPatchII のご紹介/CiPA hERG 薬物動態アッセイの QPatch での実施について ソフィオンバイオサイエンス株式会社 鶴留 一也
14.40-15.20	コーヒーブレイク
15.20-16.00	イメージングによる心筋細胞収縮評価手法-SI8000 Cell Motion Imaging System- ソニーイメージングプロダクツ&ソリューションズ株式会社 早川 智広
16.00-16.40	High-resolution <i>in vitro</i> electrophysiology system for detailed characterization of neuronal cells and networks Dr. Marie Engelene J.Obien, VP Marketing and Sales, Maxwell Biosystems AG, Switzerland
16.40-17.00	ご挨拶 シミックファーマサイエンス株式会社 代表取締役社長 小作 寛
18.00-20.00	レセプションパーティー