

Research Assistant, Japan

Do you want to be part of a growing team and play an important role in continuing our success in building the Japanese market?

Sophion was founded in year 2000 by a group of passionate electrophysiologists, all having the shared goal of making patch clamping objective and independent of user skills to provide faster, more accurate and objective results. We are an innovative, rapidly expanding company that has developed from a startup into a global organization, while maintaining the passion and drive to provide solutions for high performance cell analysis.

Today the company's analytical platforms QPatch and Qube are placed at leading pharmaceutical companies and Contract Research Organizations. With our dedicated support and service, this ensures us a leading position in the market. We try our best to be a great company to work for; fun, fulfilling and always challenging as we compete in a fast-moving market and industry.

About the position:

Due to our continued growth, Sophion Bioscience K.K. is seeking a Research Assistant to be based in our facility in Japan (Honjo Waseda, Saitama). As a Research Assistant at Sophion, you will encounter interesting challenges of technical and application character of nearly all known ion-channels. You will be challenged with the very fundamentals of patch-clamping as well as the implementation of cell cultures for various ion expressing cells and the execution of experiments using those cultured cells. You will be a part of a global application team with colleagues in Denmark, United States and China.

Main Responsibilities:

- Take care of all ion channel expressing cell lines and provide the cultured cells for patch-clamp experiments conducted on automated patch-clamp systems.
- Execute contract research experiments independently following test protocols submitted by Application Scientists and provide the scientists with test results in a timely manner.
- Maintain laboratories good conditions to make it possible to prepare chemicals, cells, consumables and solutions in a harmonious way.
- Support pre-sales support with demos and customer presentations.
- Perform quality control experiments for systems and consumables.
- Participate in validation studies for new products.
- Conduct experiments for optimizing cell lines for automated patch-clamp systems.
- Obtain test results for writing application reports, posters and publications.
- Travel as needed for customer support, user meetings, and relevant scientific conferences.

Key Qualifications:

- Experiences with mammalian cell culture and experiments using cultured cells are essential.
- Required: graduated from a university or a professional training college in science.
- Manual or automated patch-clamp experience is highly appreciated.
- Knowledge of drug discovery process or experience of in vitro contract research is valuable.
- Experience of logistic laboratory management is recognized.

Personal Qualities:

- Required: already resides in or being allowed relocating to Honjo Waseda or adjacent regions.
- Ability to work independently as an experimenter who can complete routine experiments by oneself.
- Service minded, enthusiastic, self-driven, and enjoy solving problems with a "can do, will do" mindset.
- Highly driven and motivated in work ethic.
- A hands-on person who thrives in a laboratory environment.
- Able to communicate in English with colleagues outside Japan through emails.
- Proficient in Excel, PowerPoint, Word.

Experience:

- Cell Culture (cell lines and primary cells, stem cells is a plus): 2 years' experience (preferred)
- Operation of manual and/or automated patch-clamp: 2 years' experience (ideal)
- Customer facing at CRO or drug discovery in pharma company: 2 years' experience (ideal)

Send your application and resume to job@sophion.com with "Research Assistant EMPD29251" in the subject heading.

If you have any questions, please contact the country manager in Japan:

Yuji Tsurubuchi
Country Manager, Japan
Sophion Bioscience K.K.
yts@sophion.com
Phone +81 90 2454 7967

Application deadline: As soon as possible