

Software Developer

Do you want to be part of a growing team and play an important role in continuing our success as the leading provider of automated screening equipment for preclinical drug discovery and disease modeling?

Job description:

As a software developer, you will work with software for controlling and analyzing a new system in development that executes cutting-edge biological experiments. You will join a team of scientists and engineers from different disciplines that work together to solve interesting technical and application challenges. You will be one of two software engineers that develop the software for a new product that we are currently developing. You have excellent programming skills in C++ and are adept at interfacing hardware communication protocols, databases, and programming in Windows. Familiarity with event handling and databases (SQL) is a plus.

Your integrity and ethics are high, and you have an immense focus on developing code following best practices to the highest standards. The most important skills to us are that you have a logical mind, the curiosity to learn new practices, and an ability to focus on the development of a new product within a small team.

Your key tasks:

- Develop software to control hardware, acquire and analyze data from biological experiments for a new product.
- Contribute to the design and implementation of new features for our product.
- Contribute to the definition of the global software architecture of our product.
- Participate in code reviews.
- Participate in good practices such as unit testing, automatic testing, CI pipelines.
- Participate in the product development and features definition for the following versions of our products.

Your key qualifications:

- Bachelor's degree or higher in computer science or similar field.
- 3+ years of experience in Software development in C++
- Object-oriented programming and its principles.
- Windows software development.
- Knowledge of hardware communication protocols.
- It is a plus if your profile also matches one or more of the following qualifications:
 - Hands-on experience in databases.
 - Hands-on experience in GUI development.
 - Experience with Java and C#
 - Experience with diverse version control systems.
 - Experience with JIRA and agile methodologies.
- Good communication skills in English

Sophion Bioscience A/S

Baltorpvej 154,
2750 Ballerup
Denmark
T +45 44 60 88 00
Sophion.com
info@sophion.com

Sophion Bioscience, Inc.

213 Burlington Road, Suite 105,
Bedford, MA 01730
United States
T +1 877 898 9010

Sophion Bioscience K. K.

Incubation on Campus Honjo Waseda #V203,
1011 Nishitomoda, Honjo-shi,
Saitama 367-0035, Japan
T +81 495 71 8811

Sophion Bioscience (Shanghai) Co., Ltd.

Room 550, Floor 5, Building 21
No. 1-28, Lane 588 Tianxiong Road, Pudong District
Shanghai 201318, China
T +86 21 52037529

**Your personal qualities:**

You are curious, enthusiastic and enjoy solving complex problems, and have a general technical flair. You step in when colleagues need help, and it comes naturally for you to ask for help when needed.

About us:

Sophion Bioscience is a leading global life science company founded in 2000. We are headquartered in Denmark and have subsidiaries in the United States, Japan, and China.

We specialize in developing and manufacturing automated patch clamping and cell line solutions. With our complete technical, biological, and application support, we help our partners pioneer ion channel research and drug discovery. 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 biological cell analysis.

Today, we are growing and expanding our product portfolio into emerging scientific areas that will make drugs safer for people and reduce the use of animals in drug discovery.

Sophion Bioscience is an equal-opportunity employer. We evaluate qualified applicants without regard to race, color, national origin, religion, gender, age, marital status, disability, sexual orientation, gender identity, or any other characteristic recognized by law. We do our best to be a great company to work for; fun, fulfilling, and always challenging as we compete in fast-moving markets and industries.

Please apply at:

If you think this position would be a great opportunity for you, and that you would be a great addition to the team, even though you fulfill only some of our requirements, please send your application and resume to Job@sophion.com with [Software Developer JD39427] in the subject heading.

Hiring manager:

If you have any questions, please contact hiring manager:

Sandra Wilson, Innovation & New Technology Development Manager, +45 40207753, swi@sophion.com

Application deadline:

March 25th, 2024