



## Software Engineer

### Who is Axion BioSystems?

Imagine if the scientific questions medical professionals ask were unencumbered by overcomplicated instrumentation. At Axion, we build flexible products that keep pace with today's research while striving to test the limits of one's imagination. Axion BioSystems initially launched the first commercial multi-well, Micro Electrode Array (MEA) system, the Maestro™ in 2012. Over the past 10 years, we have developed a steady pipeline of innovative, industry-leading products that increase throughput, functionality, and information for industrial and academic research. Our products are actively used to research Epilepsy, Parkinsons, ALS, toxicology, cardiac safety, and cancer with new discoveries being made constantly by our customers. If you want to be a part of a company that has the potential to save lives and shape the medical device industry, Axion may be the place for you.

### What You Would Bring to Axion BioSystems

As a member of the Axion Biosystems software team, you will contribute to the mission critical software that runs both on our physical devices and our desktop applications. Responsibilities of this role are to generate designs, author code, and review the code of other team members. You will be expected to help bring new features and products from the initial ideas to production software solutions that are used daily by researchers. To this end, you'll also work directly with Axion's multidisciplinary team to help gather, author, and maintain requirements documents for Axion's software products. As an Axion Software Engineer, you will work collaboratively with our software team, using agile methodologies with cutting-edge tools. You will be encouraged to both give and receive feedback in a supportive and constructive manner to make the team and the products we produce the best they can be.

### What We Can Offer

- Our office is in a walkable area in between Buckhead and Midtown
- We offer flexible working hours
- We have a generous benefits package including paid time off, a 401k match and bonus opportunities throughout the year
- This role reports directly to our VP of Software
- The chance to work with really smart people and cutting-edge technology solving important, life-changing problems.

### Required Skills

- Proficiency in Object-Oriented (OO) programming concepts
- Experience working in both Linux and Windows Operating Systems
- Knowledge of GIT or similar version control systems
- Experience with bug/issue tracking software
- Strong communication skills, both written and verbal





### Preferred Skills

- **Experience and Proficiency with .NET / C# / WPF**
- **Hardware Skills; including the use of oscilloscopes/signal generators, debugging connections, etc..**

