

Software Developer in Test

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 team, you will contribute to our best-in-class hardware and software. This role is responsible for the overall quality of our software products which includes firmware, desktop applications and the web. You will be writing and executing both manual and automated tests to verify the behavior and accuracy of our systems.

Our products are designed, built and assembled in our Atlanta office. 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

- A Midtown office in a walkable area
- Flexible working hours
- A generous benefits package including paid time off, a 401k match and bonus opportunities throughout the year
- The chance to work with really smart people and cutting-edge technology solving important, life-changing problems.

Duties and Responsibilities

- Write and execute tests to verify new functionality as it's developed



- Automate tests where possible so they can run automatically on any new code changes
- Follow our Software Development Process and make sure issues and bugs are handled appropriately

Required Skills

- Care deeply about the user's experience with our products and software
- Strong attention to detail
- Programming experience – ideally in C# but other object-oriented languages will do
- Familiar with Bug Tracking systems (e.g., JIRA)
- Experience working in both Linux and Windows Operating Systems
- Familiar with GIT or similar version control systems
- Strong communication skills, both written and oral

Preferred Skills

- Experience with .NET / C# programming
- Experience writing automated tests and/or automated test verification/validation

