

SENIOR SCIENTIST

What is Axion BioSystems?

Axion BioSystems, Inc. is a life science company located in Atlanta, GA. At Axion BioSystems, 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 microelectrode 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, Parkinson's disease, ALS, neurotoxicology, cardiac safety pharmacology, 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 upcoming translational therapies, Axion may be the place for you.

What would you bring to Axion BioSystems?

We are currently seeking an energetic scientist to join Axion's Applications team. Our team identifies, develops, and implements advanced applications for the Maestro family of products. Our goal is to establish and refine assays that are scientifically rigorous, yet also accessible and scale-able, including the creation of "seizures-in-a-dish" or assays for cancer immunotherapy. Specifically, this role will lead development of new analytical algorithms for existing assays in neuroscience and immunotherapy, and contribute to technical research and development for new products and applications. As a Senior Scientist on the Applications team, you would own significant development efforts from inception to delivery and routinely participate in cross-team collaboration.

Required Skills:

- PhD in Biomedical Engineering, or similar field, with 1-2 years of relevant experience, or a Master's degree with 3-5 years of relevant experience
- Experience with digital signal processing and/or computational analysis
- Proficient with one or more software languages (e.g., Matlab, Python, C#, R)
- Capable of communicating in a cross-functional engineering team
- U.S Citizenship or permanent residency required

Preferred Skills:

The following are not required, but will enhance an applicant's candidacy:

- Specialty knowledge in computational neuroscience, ranging from action potential models to neural network models
- Specialty knowledge in impedance spectroscopy
- Previous experience working with large data sets

Expectations:

The following list describes various responsibilities of this role:

- Contributes to applications development at the task, project, and team levels
- Participates in project definition, scope, planning, and execution
- Owns significant development efforts from inception to delivery
- Manages time well across multiple projects in parallel
- Able to work independently on most tasks
- Effective cross-team communication to drive projects forward

CONTACT INFORMATION:

If you are interested in this position, please send your resume and cover letter to careers@axionbio.com