



SENIOR APPLICATIONS 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. As a member of the Applications team, you would develop and maintain cell-based models, participate in the design and execution of experiments, analyze data using Axion's software tools, and produce customer-facing applications documents.

Required Skills:

- PhD in Biology, Biomedical Engineering, or similar field, or a Masters degree with 1-2 years of relevant experience
- Experience with cell-based assays in the field of immuno-oncology
- Biostatistical and data analysis experience.
- Attention to detail with good documentation and communication skills
- U.S Citizenship or permanent residency required

Preferred Skills:

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

- Specialty knowledge in the areas of genetics or high throughput screening.
- Peer-reviewed publications in the field of immuno-oncology

CONTACT INFORMATION:

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

