

## FPGA Developer

### 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. Responsibilities of this role are to design, develop and maintain FPGA firmware. You will be expected to help bring new products from the initial ideas to a finished product by working directly with Axion's multidisciplinary team. 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.

Additionally, we recognize that many FPGA developers have one or more additional skills. If your skill is hardware, you'll help with design, test and optimization of our hardware. If your skill is software, you'll work throughout the entire software development lifecycle and specify, design and implement firmware and software to run on our custom hardware. If you have other skills, let us know!

### 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.

### Required Skills

- Experience with FPGA design and the ability to be the local FPGA design expert (this means you would inform all FPGA-related decisions at Axion)
- Proficiency in the design, development, testing, debugging and maintenance of FPGA firmware in VHDL. At Axion this includes:
  - Custom peripherals to interface with high-speed ADCs and DACs
  - Digital Signal Processing and filtering
  - Ethernet Interfaces from FPGA logic
  - Embedded hard- and soft- processor SoC designs
- Experience working in both Linux and Windows Operating Systems
- Knowledge of GIT or similar version control systems
- Strong communication skills, both written and oral

### Preferred Skills

- Experience and Proficiency with .NET / C#, C or C++
- Hardware Skills; including the use of oscilloscopes/signal generators, debugging connections, etc..)
- Experience reading and reviewing board schematics
- Experience with bug/issue tracking software

