

PRODUCT DESIGN ENGINEER

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.

Axion BioSystems is seeking a Product Design Engineer reporting to the Product Design Director to design the company's next generation of products. The role will focus on designing and integrating mechanical and electrical assemblies to develop prototypes and eventually a finished product. Day to day tasks will include CAD design and review, prototype testing, accelerated life testing, process development, and documentation. Experience with low to mid volume manufacturing methodologies and best practices are critical to the role. The Product Engineer will work collaboratively across a cross functional team of engineers and scientists to solve critical problems as they relate to launching a new product.

Job Duties

- Design, build, and characterize electro-mechanical devices and subassemblies.
- Create and perform functionality testing as well as accelerated life testing on components and assemblies.
- Collaborate with other engineers and end users to work toward the best product solutions.
- Work with packaging vendors to develop custom packaging solutions to ensure safe delivery of end products.
- Create standard operating procedures, bill of materials, and build processes in preparation for manufacturing transfer.
- Create incoming inspection criteria and testing protocols for critical components.
- Develop and maintain relationships with suppliers during prototype and manufacturing transfer phases.
- Interface with consultants as needed for external expertise.
- Support established products by developing improved testing procedures and working to improve product quality.
- Attention to detail and organization skills are critical to success.

Job Requirements

- Bachelors or higher in Mechanical Engineering or related degree required.
- 2-4 years experience in an engineering or design related role.
- Experience with 3D CAD programs (CREO preferred).
- Experience generating 2D CAD drawings for fabrication.
- Understands typical manufacturing processes used to customize plastics, metals, and electronic circuit assemblies.
- Experience working in a regulated industry.

Please submit resumes to careers@axionbio.com for consideration.

