



Hardware Sustaining Engineer

Axion BioSystems, a life science company located in Atlanta, GA, is a leading manufacturer of microelectrode array (MEA) and impedance systems used to evaluate the biology of networked cells (e.g. neurons, cardiac cells, lung and liver) for applications such as cancer research, vaccine development, modeling neurological disease and drug development. When combined with modern cell culture technologies, Axion's platform allows scientists to model human biology in a dish. Axion's products are used by pharmaceutical and biotech companies, governmental agencies, and universities for drug safety, drug discovery, cell and gene therapy development and disease modeling assays.

Axion BioSystems, Inc. is seeking a Hardware Sustaining Engineer to analyze and improve upon the company's flagship Maestro multi-mode hardware platform. The ideal candidate will work well in a cross functional team where attention to detail is critical. Product improvements may include design changes to custom circuit boards, cable assemblies, custom mechanical assemblies, and the integration of off-the-shelf components. Communication with both internal and external resources to understand the impact of the changes will be important. A strong electrical engineering background with a focus on hands-on problem solving is critical. Moderate mechanical aptitude is desired.

Job Duties

- Perform root cause analysis of electronic system and component failures during manufacturing.
- Evaluate support logs to determine field failures and identify mitigating solutions.
- Improve designs by integrating test points (board level, process level, at supplier, etc) and designing out discovered failure modes.
- Identify, design, test, and integrate continuous improvement projects into manufacturing process.
- Test new processes independently with the goal of presenting improvements to Axion's Change Review Board.
- Work with suppliers to improve quality based upon processing failure analysis.
- Write and modify documentation to be followed by manufacturing technicians.
- Use Enterprise Resource Planning (ERP) tool to evaluate work in process and establish yield metrics.

Job Requirements

- BS in Electrical or Computer Engineering, or related engineering degree required.
- 2-4 years experience in Quality, Development, Sustaining, or Manufacturing related role.
- Experience with electrical test equipment (oscilloscope, multimeter), soldering, electronic debugging.
- Ability to communicate problems at both an executive and detailed level.
- PCB Layout and Schematic Capture software experience.
- QMS experience and proper documentation practices desired.
- Self-motivated.
- Work is primarily performed on-site in Axion's Atlanta office.

Excellent references and a demonstrated record of success are essential. A resumé along with a letter of introduction should be sent to careers@axionbio.com.

