

## Senior Director, Operations

Axion BioSystems, is a Private Equity backed life science tools company located in Atlanta, GA. The Company 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 Senior Director, Operations reporting to the CEO and responsible for managing, coordinating, and improving the company's Operations team. The ideal candidate will have diverse experience managing all aspects of a successful operations team in a rapidly growing company including supply chain, site identification, production, shipping/receiving, and warehousing/inventory management. Excellent problem solving skills are critical to the position to resolve challenges as they arise. Effective personnel management is also essential to successfully manage this growing team.

### Job Duties

- Direct and manage all Operations team functions and personnel.
- Organize production plans and work on improving output efficiency.
- Assist with supply chain forecasting and decision making as issues arise, including the identification of new suppliers and cost saving opportunities with remote suppliers.
- Help to identify new production site and plan layout and move of the operation.
- Work with Manufacturing Engineering team to identify and implement opportunities to improve productivity.
- Develop relationships and develop contracts with key suppliers.
- Set and execute on goals, key metrics, and output targets for the team.
- Coordinate with Customer Success team to schedule customer system shipments and inform schedule changes.
- Work with Sales to organize in-house and in-field repairs and preventative maintenance.
- Maintain safety standards and routine maintenance of company facilities.
- Utilize ERP and MES systems to track inventory, perform counts, and organize production efforts.

### Job Requirements

- Bachelors or higher in Engineering, Management, or related degree required.
- 10-15 years experience in an operations or production management related role.
- Experience with ERP and MES systems.
- Experience creating and implementing documentation, quality and production systems.
- Experience managing a diverse team.
- Excellent problem solving, verbal and written communication skills.

Excellent references and a demonstrated record of success are essential. A resumé along with a letter of introduction may be sent to [careers@axionbio.com](mailto:careers@axionbio.com).

