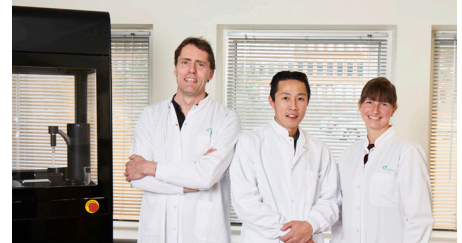


Production Engineer (PTA) - Copenhagen

Do you want to be part of a growing team and play an important role in continuing our success?

Job description:

As our new production engineer (PTA), you will be responsible for ensuring and continuously improving our high-quality level. Furthermore, with direct reference to the Head of Supply Chain and Manufacturing, you will be responsible for several key activities within our productions' technical area e.c. setting up serial production, creating uniform production documentation etc.



You have an energetic personality with strong electric or electromechanical skills. Your integrity is high, and you have an eye for quality. Experience with the production of complex electromechanical products is a plus but not a must. Being a hands-on person is your most important skill to us, and you are not afraid of taking the initiative when needed. To be a team player in our organization is a must!

Key Qualifications

- Required: Electric or Electromechanical education on bachelor level or equivalent
- Background: 3-5 years of experience with the production of smaller series of complex electrical and electromechanical products.
- Experience with analyzing RMA item to improve quality.
- Experience with functional testing of electrical and mechanical production is required
- Solid understanding of electronic design and production.
- Knowledge about structured DFM on electrical and mechanical modules.

Main Responsibilities

- Managing production and production processes of electromechanical instruments.
- Bridging products released from R&D to production
- Plan and perform incoming and outgoing quality control
- Setup test procedure for PCB, IPC, modules etc
- Write and maintain working instruction and guides.
- Ensure quality measures are anchored inside the production

Personal Qualities

- Ability to work independently and prioritize in a dynamic working environment
- Service-minded, enthusiastic, self-driven, and enjoy solving complicated problems with a general technical flair and a "can do, will do" mindset.
- Able to communicate fluently in English. Understanding Danish is an advantage.
- Able to step in when colleagues need help, and not afraid to ask for help when needed
- A hands-on person who thrives in a development environment
- Proficient in Excel, PowerPoint, and Word

Send your application and resumé to job@sophion.com with the job title "Production Engineer" in the subject heading.

If you have any questions, please contact the recruiting manager:

Jesper Haslund
jha@sophion.com
Phone +45 30 73 84 47

Application deadline: June 13th, 2021

Sophion was founded almost twenty years ago by a group of passionate electrophysiologists, all having the shared purpose of making patch clamping objective and independent of user skills to provide faster, more accurate and objective results. We are an innovative, rapidly expanding company that has developed from a startup into a global organization, while maintaining the passion and drive to provide solutions for high-performance cell analysis.

Today the company's analytical platforms QPatch and Qube are placed at leading pharmaceutical companies and Contract Research Organizations (CRO) and with our dedicated support and service, it ensures us a leading position in the market. We try our best to be a great company to work for; fun, fulfilling and always challenging as we compete in a fast-moving market and industry.