



ARBOR ASSAYS
Interactive Assay Solutions™

Are you an organized and detail-oriented individual who has **2+ years of experience in a production/manufacturing environment?** Can you troubleshoot equipment and recommend innovative process improvements? Do you want to work for an established and respected company, the **first employee-owned life sciences company in the world** that will provide you with opportunities for professional growth? If so, please read on!

A DAY IN THE LIFE OF A MANUFACTURING ASSISTANT

As a Manufacturing Assistant, you work as part of a team in a lab environment to support various manufacturing functions to include filling, capping, labeling and packaging products in support of the production schedule.

With attention to detail, accuracy, and timeliness, you:

- **Ensure pertinent documentation is in-place and filled out.**
- **Carry out and maintain inventory control measures.**
- **Spend time participating in inventory verifications and cycle counting.**
- **Provide back-up support to our Shipping Coordinator.**

WHAT YOU NEED TO SUCCEED AS A MANUFACTURING ASSISTANT

- Minimum of high school diploma and 2 years of experience in a production environment. Experience in life sciences industry preferred.
- Detail-oriented, highly responsible and possess excellent organizational and communication skills.
- Ability to troubleshoot equipment and recommend innovative process improvements.
- Ability and desire to follow standard procedures without deviation.
- Ability to work independently and collaboratively in a production team.
- Represent the company's best interests, mission and community focus.
- Must be able to perform occasional lifting up to 30 lbs. individually and 50 lbs. with assistance.

The Manufacturing Assistant position earns a **competitive wage**. We also offer **medical, dental and vision effective your first day**. Along with an IRA and voluntary life insurance. If this sounds like the right opportunity for you, apply today!

ABOUT ARBOR ASSAYS

We are a life sciences company based in Ann Arbor, MI. Founded in 2007 and employee-owned since 2017, Arbor Assays has always had the simple desire to create the highest quality assay products for the research community, using US-sourced materials and a pro-environment perspective. We take great pride in building products that make someone's job easier, help create sound research and provide consistent results.

BE A PART OF A COMPANY THAT IS GOING PLACES. JOIN ARBOR ASSAYS!