



ARBOR ASSAYS
Interactive Assay Solutions™

Are you a Research Lab Technician who has 1+ years of experience with ELISA or PCR assays? 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!

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. Alongside our enthusiasm for science and research, **we have a passion to see each and every one of our employees grow and succeed both personally and professionally.**

A DAY IN THE LIFE OF A RESEARCH LAB TECHNICIAN

As a Research Lab Technician, you work as part of a team in a lab environment to complete client-driven projects. With attention to detail, accuracy and timeliness, you assist in ELISA, EIA, Detection assay development experiments using standard biochemistry techniques. You also assist in protein purifications, protein/ small molecule extraction, UV/Vis analysis and standard reagent preparation. Following established SOPs, you perform sample analyses. This includes maintaining a detailed and accurate record of your experiments and updating R&D lab notebooks, SOPs and general R&D documentation. As a Research Lab Technician, you will also spend some of your time maintaining a general consumable lab supply and assisting with Quality Control and Manufacturing activities.

WHAT YOU NEED TO SUCCEED AS A RESEARCH LAB TECHNICIAN

- Bachelor's degree and 1+ years of experience in the life sciences industry.
- Previous experience with biological sample handling at a lab highly desired.
- Detail-oriented, highly responsible, and possess excellent organizational and communication skills.
- Ability and desire to follow standard procedures without deviation.
- Ability to apply critical thinking for suggested efficiency changes.
- Ability to work independently and collaboratively in a research team.
- Desire and ability to collaborate to meet company goals and objectives.

This Research Lab Technician position earns a **competitive salary**. We also offer a **full suite of benefits** to choose from, including **medical, dental, and vision effective your first day. A 401(k) plan, and voluntary life/ADD&D insurance.** If this sounds like the right opportunity for you, apply today!

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