



**Title:** QC Specialist – Biochemical Assays

**Position Description:**

Reaction Biology (RB), is a private pharmaceutical research company. We provide contract research services to over 1500 clients worldwide, through our labs in Malvern, PA and Freiburg, Germany. We are currently seeking a QC Specialist for our Malvern, PA facility. We are looking for a reliable and motivated individual with a strong science background to work in our lab.

The successful candidate will provide quality control review of laboratory records, reports, and other documentation for accuracy, appropriateness, and timeliness.

**Job Responsibilities:**

- QC data in an accurate and timely manner, balancing multiple service orders, at times in high volumes.
- Work with lab scientists to meet customer demands while always meeting established quality control standards.
- Provide knowledgeable and accurate feedback regarding data quality to Research Associates and Research Scientists.
- Support Research Scientists and Research Associates and help them as required, including assisting in troubleshooting issues that arise in the data.
- Help to provide oversight over the QC process.
- Work with lab leadership to develop, revise, implement, and maintain standard operating procedures (SOPs)
- Follow SOPs and maintain good documentation during work activities.
- Working in a team environment to ensure right on time first time delivery for clients.
- Performing training of new staff and maintains training documentation. Re-assessing and re-training staff found to be involved in steps that resulted in errors.
- Organizing and participating in calls and virtual meetings with scientists and managers as needed.

**Required Knowledge, Skills, and Abilities:**

- Bachelor's degree (required) in a related field with 2+ years of experience or Master's degree in a related field with 1+ years of experience.
- Experience working in QC/QA preferred.
- Previous experience with enzymatic assays is desired
- Attention to detail, and organization skills are critical.

Let's discover together.



- Collaborative and proactive attitude with excellent written and oral communication skills.
- Flexibility to accommodate to rapidly changing priorities and deadlines in a fast-paced and dynamic team-based environment.

The work environment is casual; however, professional behavior, a collaborative attitude, attention to detail, and the ability to meet deadlines are absolute requirements.

Finally, all employees are obligated to meet the highest standards with respect to honesty, accuracy, conflict of interest, intellectual property rights, confidentiality and respect for information sources and readers.

Benefits include health insurance, 401-K plan, and paid vacations. RB is an equal opportunity employer.

