

Research Associate II – In Vitro Job Posting

Position Summary:

The In Vitro Research Associate II is responsible for conducting cell culture and in vitro studies to support drug development and research under general supervision. This role is responsible for performing laboratory tasks with precision while adhering to established protocols, regulatory guidelines, and study objectives.

Main Responsibilities:

- Conduct in vitro experiments and assays following established protocols and biosafety guidelines.
- Prepare reagents, cell cultures, and other necessary materials for experiments.
- Assist in the development, optimization, and validation of in vitro models and assays.
- Operate independently with minimal supervision on varied and occasionally difficult assignments, adhering to standard operating procedures (SOPs) and study-specific documents (SSDs) to ensure compliance and accuracy in study execution.
- Take full ownership of assigned studies/assays, independently managing all aspects of execution, from planning through completion, without needing constant direction or oversight.
- Set priorities and organize work within general guidelines established by management and in consultation with more senior technical staff.
- Independently identify and prepare the necessary supplies and materials required to complete assigned tasks and studies.
- Operate independently with minimal supervision on varied and occasionally difficult assignments, adhering to standard operating procedures (SOPs) and study-specific documents (SSDs) to ensure compliance and accuracy in study execution.
- Take full ownership of assigned studies/assays, independently managing all aspects of execution, from planning through completion, without needing constant direction or oversight.
- Set priorities and organize work within general guidelines established by management and in consultation with more senior technical staff.
- Assist in the preparation and maintenance of study documentation to ensure compliance with protocols and regulatory requirements.
- Contribute to the Quality System by participating in audits and supporting the preparation and review of SOPs.
- Independently identify and prepare the necessary supplies and materials required to complete assigned tasks and studies.
- Other responsibilities/tasks as assigned.

Key Qualifications:

- Bachelor's degree in biology, biochemistry, or a related scientific discipline.



- Minimum of 3 years' industry experience.
- Proficiency in Microsoft Office, including Office 365 applications and data entry, is preferred.
- Strong attention to detail with a commitment to accuracy and precision.
- Excellent interpersonal and communication skills.
- Proven ability to work effectively in a team-oriented, collaborative environment.
- Previous industry experience preferred.

Physical Demands:

- Frequent standing, walking, bending, and reaching to perform animal handling, dosing, and sample collection.
- Lifting, carrying, and moving equipment and supplies.
- Dexterity and fine motor skills required.
- Prolonged periods of wearing personal protective equipment (PPE).
- Ability to work flexible schedules, including weekends and holidays.
- Visual acuity and attention to detail.

Reaction Biology Corporation is an Equal Opportunity Employer that does not discriminate on the basis of actual or perceived person's race, color, national origin, age, religion, disability status, sex, sexual orientation, gender identity or expression, genetic information or marital status, or any other characteristic protected by applicable federal, state or local laws. Reaction Biology is an E-Verify employer.