

## **Research Associate II – In Vivo Job Posting**

### **Position Summary:**

The In Vivo Research Associate II is responsible for conducting in vivo studies to support drug development and research under minimal supervision. This role ensures the proper care and welfare of research animals while following protocols, regulatory guidelines, and study objectives.

### **Main Responsibilities:**

- 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.
- Conduct in vivo studies, including tumor implantation, surgical procedures, compound administration, and biological sample collection, while accurately recording and reporting data.
- Perform luminescent and fluorescent imaging to assess drug distribution and monitor disease progression in animal models.
- Conduct downstream analyses of ex vivo samples to support study objectives.
- Assist in the preparation and maintenance of study documentation to ensure compliance with protocols and regulatory requirements.

### **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.