



Job Title	Research Associate I	Reports to	Department Manager
Department	Biochemical	Location	Malvern
	Non-Exempt	Date	16 January 2024

Reaction Biology is a global contract research organization (CRO) with laboratory facilities in Malvern, PA, Hershey, PA, Freiburg, Germany and Heidelberg, Germany. We provide drug discovery research and development services to biopharmaceutical customers worldwide. Our capabilities and services include protein & peptide production, biochemical and biophysical assays, high-throughput screening, an extensive array of cell-based and pharmacological assays modeling various diseases, exploratory toxicology, biomarker discovery, and pharmaceutical batch release testing.

Since its inception in 2001, Reaction Biology has grown to become a global leader in drug research and development, assisting in the discovery of life-saving therapeutics for diseases such as cancer and neurodegeneration. Reaction Biology has assisted in the advancement of many of its customers' drug candidates through the preclinical pipeline stages into later clinical phases. Reaction has one of the world's largest panels of kinase assays, with over 780 kinases, and performs more than 5,000 client projects annually with over 2,000 validated assays and growing.

#### **Position Summary:**

To provide ground level lab work following standard protocols according to department needs. The employee will support other team members in the execution of routine lab procedures, as well as to help sustain the functionality of the lab through regular cleaning, stocking inventory, and maintenance.

#### **Main Responsibilities:**

- Prepare and handle reagents.
- Maintain radiation safety.
- Maintain lab inventory.
- Safely handle client enzymes.
- Basic plate reader operation.
- Customer source plate preparation.
- Maintain lab safety – follow proper waste disposal.
- Maintain electronic logs.
- Basic lab equipment use, as determined by department.
- Understanding and use of Reaction Biology specific programs for sample, data, and order tracking
- Perform standard lab calculations.
- Use proper pipetting techniques.
- Perform data analysis, as expected by department.
- Perform proper lab and office maintenance and cleaning – practice good lab practices, follow daily and weekly cleaning checklists.
- Clearly communicate verbally and in writing.
- Other duties as assigned.



**Key Qualifications:**

- Bachelor of science degree.
- Proficiency in lab setting.
- Basic understanding of lab equipment
- Ability to work independently and in teams.

**Physical Demands:**

- Sedentary work that primarily involves sitting/standing
- Light work that includes moving objects up to 30 pounds

Reaction Biology Corporation is an Equal Opportunity Employer that does not discriminate on the basis of actual or perceived person's race (including hairstyle/texture), 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.