



Title: Research Associate III – Biophysical Assays

Job 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 looking for a reliable and motivated individual with a strong science background and experience with SPR analysis to join a team involved in early stage drug discovery studies. The position is a bench position with primary duties in executing a variety of biophysical assays with a focus on SPR analysis.

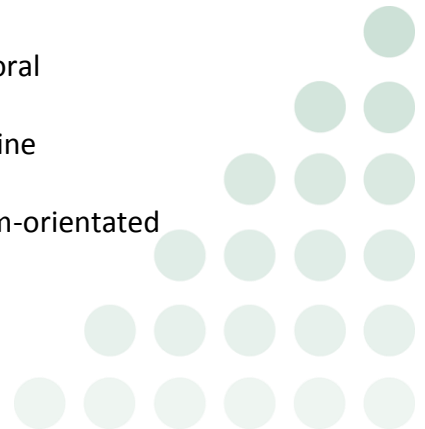
Job Duties and Responsibilities:

- Maintain lab inventory for the biophysical assay service group
- Complete various SPR assays for external customers with high-quality data and in a time-efficient manner
- Establish new SPR assay platforms
- Manage multiple projects and deliver timely results
- Screen and characterize compounds or other type of molecules and determine the mode of molecular action
- Be able to utilize literature and internal data sets to design and test target modulation hypotheses
- Assist with administrative duties of sending/receiving client forms
- Help maintain a calendar system for scheduling of projects and deliverables

Required Knowledge, Skills, and Abilities:

- BS or equivalent in biochemistry, biophysics or related discipline, with 1-2 years of wet lab experience in a relevant discipline
- MS or equivalent in biochemistry, biophysics or related discipline, with 0-1 year of wet lab experience in a relevant discipline
- Detail-oriented
- Collaborative and proactive attitude with excellent written and oral communication skills
- Flexibility to accommodate rapidly changing priorities and deadline
- Ability to work in a team-based environment
- Ability to work independently with minimal supervision in a team-orientated environment

Let's discover together.



Preferred Knowledge, Skills, and Abilities:

- Hands on experience with SPR and/or other biophysical tools (MST, ITC, NMR, etc.)
- Proven expertise in assay development using biophysical techniques
- Familiarity with SPR theory and kinetic analysis
- Experience in kinetic profiling and library screening of small molecules is a plus

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 at RB include health insurance, 401-K plan, student loan benefit, and paid vacations. RB is an equal opportunity employer.

