



Title: Senior Scientist – 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 seeking 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.

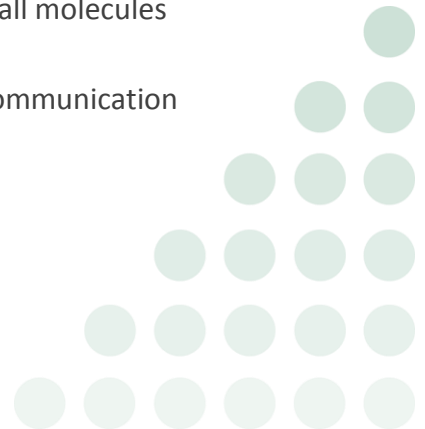
Job Duties and Responsibilities:

- 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
- Interact directly with customers in order to understand their requests, to trouble shoot assay problems and to propose/pursue solutions
- Screen and characterize a variety of molecules and determine the mode of action with a primary focus on low molecular weight analytes
- Be able to utilize literature and internal data sets to design and test target modulation hypotheses

Required Knowledge, Skills, and Abilities:

- PhD or equivalent in biochemistry and biophysics or related discipline, with 2+ years of SPR working experience
- Demonstrated in-depth understanding of SPR theory and proficiency in assay development, optimization, kinetic profiling and library screening of small molecules
- Attention to detail
- 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

Let's discover together.



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

