



Title: Research Associate – 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 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 binding kinetics determined via an optical technique known as surface plasmon resonance (SPR).

Job Duties and Responsibilities:

- Maintain lab inventory for the biophysical assay service group
- Prepare buffers and other reagents for use in experiments
- Participate in cleaning procedures for lab equipment
- Help maintain the SPR instruments via weekly and monthly cleaning procedures
- Assist in the preparation of 96- and 384-well plates using a multi-channel pipette to perform serial dilutions of samples following a set-protocol
- Help with establishing SPR assays for new targets by following a general protocol to explore different buffer conditions and sample preparations
- 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

Let's discover together.



Preferred Knowledge, Skills, and Abilities:

- Experience with multi-channel pipettes
- Experience preparing a variety of buffers and lab reagents
- Experience with SPR (or other biophysical techniques) would be a plus

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

