



Title: Senior Scientist – Biochemical Assay Development

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 currently seeking a Senior Scientist for our Malvern, PA facility. We are looking for a reliable and motivated individual with a strong science background to work in our lab.

The successful candidate will independently plan and execute internal and external research service projects in an accurate, reproducible, and timely manner.

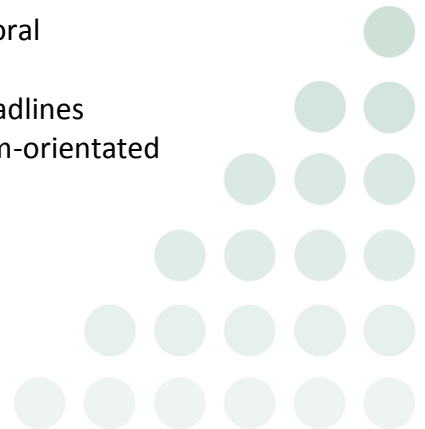
Job Duties and Responsibilities:

- Design, implement, and execute a wide variety of in vitro assays for enzyme targets.
- Screen and characterize small molecule compounds and determine the mode of compound action.
- Help to troubleshoot issues that arise in the lab and propose or pursue solutions.
- Utilize literature and internal data sets to design and test target modulation hypotheses.
- Maintain and further develop new and existing scientific relationships with our clientele by helping to provide a consultative experience.

Qualifications:

- PhD in biochemistry or related discipline with 2+ years of experience
- Experience in steady-state kinetic characterization of enzyme, substrate, and inhibitor mechanisms preferred.
- Scientific and technical background in drug discovery, with demonstrated proficiency in assay development, optimization, and screening with small molecules.
- Attention to detail, organization skills, and a high degree of scientific integrity are critical
- Collaborative and proactive attitude with excellent written and oral communication skills.
- Flexibility to accommodate to rapidly changing priorities and deadlines
- Ability to work independently with minimal supervision in a team-orientated environment

Let's discover together.



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.

