



Title: Senior Scientist – Receptor Pharmacology

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.

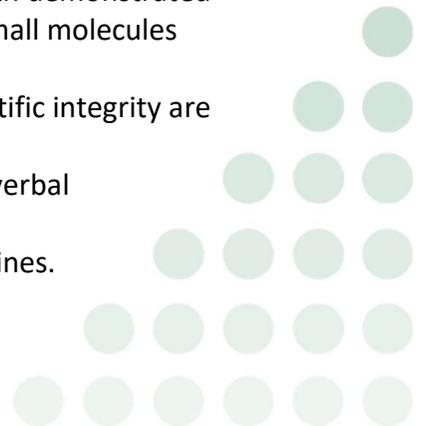
Job Duties and Responsibilities:

- Design, implement and execute a wide variety of in vitro assays for receptor (GPCR, nuclear receptor, ion channel) 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/pursue solutions.
- Utilize literature and internal data sets to design and test target modulation hypotheses.
- Maintain and further develop scientific relationships with our client base through expert consultation when needed.

Qualifications:

- PhD in pharmacology or related discipline with 0-2 years of experience.
- Experience with standard in vitro techniques including, but not limited to, ligand binding assays (kinetics, standard steady state receptor properties, SAR), standard cell signaling (calcium mobilization, cAMP modulation, beta arrestin biology) preferred.
- Preference will be given to candidates with practical experience in ion channel biology including calcium signaling and QPatch instrumentation.
- Scientific and technical background related to drug discovery with demonstrated proficiency in assay development, optimization and screening small molecules preferred.
- Attention to detail, organization skills and a high degree of scientific integrity are essential.
- Collaborative and proactive attitude with excellent written and verbal communication skills required.
- Flexibility to accommodate rapidly changing priorities and deadlines.

Let's discover together.



- Ability to work independently with minimal supervision in a team-orientated environment.

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.

