

Senior Scientist Mol and Cell Biologist

Do you want to explore frontiers of biology to make breakthrough discoveries and use its applications in drug discovery and diagnostics? Ubiquitin proteasome system (UPS) provides a challenge as a new wave of drugs are emerging, such as Protein Targeting Chimeras (PROTACs) that target undruggable space. Senior Scientists are expected to develop innovative technologies to understand the role of ubiquitin in cell physiology. Our technologies are used worldwide by academic, pharma and biotech companies. The Senior Scientist will lead research programs to develop innovative technologies, execute drug discovery projects for our clients, as well as drive the development of new products and product applications. He/she will write scientific articles, patents, and present data at national and international meetings. Please consult our website www.lifesensors.com more details.

Position:

Successful candidates will have experience in molecular and cellular biology, biochemistry molecular cloning, development of assays for UPS enzymes, protein expression, characterization and purification of proteins. Experience with UPS enzymes function is an advantage but not essential.

Required Skills, Experience and Keys to Success:

- Ph.D. in biological sciences
- 3-10+ years of laboratory experience in an academic or industrial setting
- Experience in molecular cloning and designing expression vectors
- Experience in protein expression, purification and protein engineering
- Knowledge of mass spec and biophysical techniques to study protein structure and function
- Experience in assay development and HTS
- Knowledge of the ubiquitin proteasome pathway is desirable
- Ability to work in a highly entrepreneurial environment
- Write scholarly papers, patents and provide intellectual leadership for the company
- Proven interpersonal skills and the ability to work as a team

Benefits:

LifeSensors employees enjoy an outstanding total compensation package, including excellent salaries based on qualifications and experience, and benefits including:

- Medical, dental, and vision coverage
- Generous Paid Time Off policy, in addition to standard Holidays
- 401K Retirement Savings Plan with a company match
- Bonus and Life Insurance programs
- Relocation Assistance

If you would like to play a critical role in ubiquitin research and in the drug discovery field, please e-mail your resume to careers@lifesensors.com