

## Entry Level Scientist

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. The Scientist will join a team that develops innovative technologies, execute drug discovery projects for our clients, as well as drive the development of new product applications. He/she will write scientific articles, patents, and present data at national and international meetings. Please consult our website [www.lifesensors.com](http://www.lifesensors.com) more details.

### Position:

We are seeking individuals who are passionate about research, bring out of the box thinking and take bold steps to develop innovative technologies with applications in the ubiquitin research field. Successful candidates will have basic knowledge of biochemistry and molecular biology techniques, in particular experience with protein expression and purification and cloning. Knowledge of the ubiquitin proteasome pathway is desired.

### Required Skills and Experience:

- BS or MS in biological sciences or related degree
- 0-2 years of laboratory experience in an academic or industrial setting
- Experience in biochemistry and molecular biology
- Ability to work in a highly entrepreneurial environment
- Ability to adapt quickly to change and manage multiple priorities
- 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
- Tax-effective reimbursement accounts for child care and medical expenses
- Generous PTO
- Retirement Savings Plan with a company match
- Bonus and Life Insurance programs

If you would like to play a critical role in UPS and PROTAC drugs, and product development, please e-mail your resume to [careers@lifesensors.com](mailto:careers@lifesensors.com)