

Senior Scientist Proteomics

LifeSensors has established its leadership position by providing innovative research tools for the ubiquitin pathway and SUMO-fusion protein expression platform. Scientists at LifeSensors are developing new technologies to understand the role of ubiquitin in cell physiology. Our products are used worldwide by academic, pharma and biotech companies. LifeSensors is located in the Great Valley Corporate Center, a biotech hub 35 miles west of Philadelphia, in Malvern Pennsylvania, USA.

Position:

Ubiquitylation is the next frontier in drug discovery research. To understand multitudes of ubiquitin modification of proteins mass spectroscopy applications are vital. Successful candidates will have mass spectrometry experience as well as broad experience in biochemistry and molecular biology techniques, in particular experience with ubiquitin pathway enzyme function, protein expression, purification and characterization. Experience in assay development, high throughput screening (HTS) and/or protein microarrays would be desirable.

Required Skills, Experience and Keys to Success:

- MS or Ph.D. in biological sciences
- 0-10+ years of laboratory experience in an academic or industrial setting
- Mass spectrometry experience
- Experience in biochemistry and molecular biology is required
- Knowledge of the ubiquitin proteasome pathway is desirable
- Ability to work in a highly entrepreneurial environment
- Ability to adapt quickly to change and manage multiple priorities
- 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
- Tax-effective reimbursement accounts for child care and medical expenses
- 10 days annual vacation, 10 paid holidays each year , 5 paid sick days per year
- 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