

Entry Level Scientist

LifeSensors, Inc. discovers and develops enabling technologies for drug discovery and research markets. LifeSensors has established its leadership position by providing innovative research tools for the ubiquitin pathway and SUMO-fusion protein expression platform. Our products are used worldwide by academic, pharma and biotech companies. LifeSensors is located 35 miles west of Philadelphia, in Malvern Pennsylvania, USA.

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
- 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 product production, please e-mail your resume to careers@lifesensors.com