

## General Manager, R&D

### Position Summary:

Create value by developing innovative products and technologies with applications in the Ubiquitin proteasome system (UPS) and Protein Targeting Chimeras (PROTACs) that target undruggable space. Lead a highly motivated team of junior and senior scientists in the execution of LifeSensors R&D strategic plans. Execute translation of innovative ideas into marketable technologies that increase revenue. Support marketing of LifeSensors' products and delivery of contract services with technical guidance.

### Required Skills, Experience and Keys to Success:

- Ph.D. in biological sciences
- Proven track record of 5 - 20 years in R&D at Biotech or Pharma or academic institute
- High-level understanding and experience with current biochemistry, protein chemistry and/or molecular biology technologies
- Excellent interpersonal and communication skills
- Ability to lead, strategically plan, and manage highly dynamic activities in R&D

### Essential Duties & Responsibilities:

- Provide day-to-day leadership in R&D that executes LifeSensors' strategic plans
- Manage research programs leading to novel products and technologies
- Coordinate with R and D teams to deliver milestones for internal and external projects
- Coordinate with marketing team to launch new products
- Establish collaboration with key opinion leaders to validate LifeSensors technologies
- Represent LifeSensors at national and international conferences
- Write articles in scholarly and trade journals to enhance visibility of LifeSensors and to promote our innovative technologies for UPS and PROTAC fields.

### Benefits:

A competitive compensation package including base salary that is commensurate with experience will be offered. LifeSensors offers a complete benefits package including healthcare, dental and vision insurance, 401K plan contributions, life and disability insurance, and the opportunity to work in a professionally rewarding, entrepreneurial environment that fosters personal development.

### About Lifesensors, Inc:

LifeSensors, Inc., a biotechnology company that discovers and develops enabling technologies for therapeutic, drug discovery and research markets, has established its leadership position by providing innovative research tools for UPS and PROTAC drugs. Its prokaryotic and eukaryotic protein expression platforms SUMOpro and SUMOstar are used worldwide by academic and government institutions, as well as pharmaceutical and biotechnology companies. LifeSensors is located in the Great Valley Corporate Center, a biotech hub 35 miles west of Philadelphia, in Malvern Pennsylvania, USA.

Interested candidates should submit a cover letter and resume to [careers@lifesensors.com](mailto:careers@lifesensors.com).