



SENIOR RESEARCH ASSOCIATE - GPCR ASSAY DEVELOPMENT

Montana Molecular is an established biotech company in Bozeman MT that produces live cell assays for therapeutic discovery and basic research. A core value at Montana Molecular is the belief that success depends upon maintaining a balance between scientific discovery, health, and the everyday joy of life. Our mission centers on creating advanced tools for interrogating signaling in living cells, by combining next generation genetically-encoded fluorescent biosensors with kinetic signaling analysis to produce high quality data and insight into the mechanisms of disease and cell physiology. A highly motivated cell biologist/assay developer that enjoys working in a fast-paced, results-oriented team is invited to apply for this position. Interested candidates are encouraged to submit resume and cover letter to info@montanamolecular.com.

Requirements

- 5+ years of relevant laboratory experience.
- A record of achievement in academic or industry research.
- Expertise in aseptic cell biology techniques.
- Understanding of standards for cell based assay development.
- Experience with cell signaling measurements.
- Excellent written and verbal communication skills.
- Ability to follow and implement standard operating procedures.
- Familiarity with molecular biology techniques.
- Live cell imaging and microscopy expertise is a plus.
- BS or higher degree in a relevant discipline (e.g., biology, chemistry, pharmacology).

Responsibilities

- Develop, optimize, and document GPCR kinetic signaling assays in living cells.
- Analyze and present results at team meetings.
- Prepare data for peer-reviewed scientific publication.
- Work effectively with team members to develop solutions in a customer-focused environment.
- Implement standard operating procedures.
- Maintain a detailed lab notebook on a daily basis.
- Report to the Associate Director of Assay Development and Services.

Highly competitive salary and fringe benefits package. Frequent opportunities for career development and advancement.