

FULL TIME GPCR SCIENTIST

Summary

Montana Molecular is an established biotech company in Bozeman MT. A core value at Montana Molecular is the belief that success depends upon finding the important balance between health, science, and the every day joy of life. Our mission centers on creating advanced tools for interrogating signaling in living cells. We combine next generation, genetically-encoded fluorescent biosensors with kinetic signaling analysis to produce new assays for basic research and drug discovery. A highly motivated GPCR Scientist that enjoys working on a fast-paced, results-oriented team is invited to apply.

Requirements

The ideal candidate will have:

- 3+ years of applicable laboratory experience with a record of achievement in GPCR research
- Outstanding cell biology and molecular cloning skills
- Joy in beautifully designed and executed experiments
- Expertise using software tools such as Prism, SnapGene, and Fiji
- Understanding of standards for assay development
- Excellent written and verbal communication skills

Responsibilities

- Independently optimize, and document GPCR kinetic signaling assays in living cells
- Analyze and present results at group meetings and scientific conferences
- Interact with GPCR researchers across the globe to transfer biosensor assay technologies
- Prepare data for peer-reviewed scientific publication
- Produce high quality web-based content summarizing scientific publications relevant to GPCR signaling in living cells
- Work effectively with team members to develop and implement solutions in a customer-focused environment
- Follow established lab operating procedures
- Occasionally work outside of regular business hours

Physical Demands

While performing the duties of this job, the employee is regularly required to sit for prolonged periods of time. The employee is frequently required to talk and hear; use hands to handle or feel; reach with hands and arms above shoulder height and below the waist. Using proper lifting techniques, the employee must be able to lift up to 30 lbs.

Competitive salary and fringe benefits package. Frequent opportunity for career development and advancement.

Qualified applicants may apply by sending cover letter and resume to info@montanamolecular.com