

Research Associate II, Genomics (Carlsbad CA)

Active Motif, Inc. is a privately held biotechnology company that develops, manufactures and supplies innovative cell & molecular biology-based reagents, kits and services that help researchers elucidate the function, regulation and interactions of epigenetic and nuclear processes. We aspire to provide customers with high-quality products, complemented by superior service and support.

We are seeking a highly motivated, team-orientated individual to join the Research and Development team as a Research Associate II at our Carlsbad facility. The candidate will work closely with other R&D team members to develop new genomic technologies and assays for the reagent market.

In the performance of their respective tasks and duties, all Active Motif employees are expected to conform to the following:

- Perform quality work within deadlines with or without direct supervision.
- Communicate with co-workers, customers and consumers in a courteous and professional manner.
- Work effectively as a team contributor on all assignments.
- Work independently while understanding the necessity for communicating and coordinating work efforts with other employees and organizations.
- Suggest and implement improvements to internal processes to increase productivity and streamline department workload.
- Perform other duties as required and / or assigned by management.

Responsibilities / Duties / Functions / Tasks

- Ability to plan, design and execute well-controlled experiments. Develop basic experimental protocols for routine use. Ensure availability of materials and equipment to carry out experiments.
- Make detailed observations, analyze data and interpret results. Maintain complete and thorough lab notebook. Suggest additional experiments based on research findings.
- Author draft versions of manufacturing documentation required for efficient product transfer and release to market.
- Contribute ideas and suggestions to improve standard laboratory techniques, improve protocols, processes and equipment.
- Maintain cell culture facilities and culture cell lines as required for experiments.
- Maintain current lab supplies of common reagents and re-order as needed to maintain workflow.
- Maintain familiarity with current scientific literature and competitive products.
- Perform literature reviews directly associated with assigned project(s).
- Work under immediate supervision of, and report to, a group leader.
- Other duties as required and / or assigned.

Qualifications / Requirements

- Bachelor's degree required (Masters preferred) in Biology, Biochemistry, Genomics, or other related discipline.
- Three years laboratory experience in an industrial setting with a working mastery of experimental design.
- Demonstrated proficiency in the use of basic laboratory techniques, equipment and materials and data analysis / interpretation.
- General familiarity with genomics-based assays such as chromatin immunoprecipitation (ChIP), DNA methylation assays and Next Generation Sequencing.
- Occasional domestic and international travel may be required.

Preferences

- Proficiency in the fields of epigenetics and genomics.
- Experience with ChIP, Next Gen Sequencing sample preparation.
- Experience in the analysis of whole-genome Next Gen sequencing data.

Work Requirements

- May have to work overtime / weekends to meet project deadlines.
- Ability to lift up to 50 lbs.
- May be required to sit and / or stand for long periods while performing duties.
- Must be capable of working safely with hazardous chemicals and biologicals.

In addition to a highly interactive and stimulating work environment, Active Motif also provides competitive salaries, stock options and an excellent benefits package. Active Motif is an equal opportunity employer and promotes diversity throughout its workforce. Active Motif is headquartered in Carlsbad, California, and has additional offices in Belgium, Japan and China.

If you meet the above requirements and are interested in a career with Active Motif, please submit a cover letter and resume outlining your qualifications and experience to: hr@activemotif.com, referencing "Research Associate II, Genomics 2018" in the Subject line.

We thank all applicants, however, only those selected for interviews will be contacted.