

## Research Associate II – Services (Carlsbad, CA)

Active Motif, Inc. is a privately held biotechnology company focused on supplying innovative kits and reagents for epigenetic and nuclear function research. Active Motif's Services group, which performs specialized epigenetics-based experiments for leading academic groups and pharmaceutical companies, is seeking a Research Associate to join our team.

The Research Associate will work closely with other members of the epigenetic services team at our Carlsbad facility to generate high-quality data for many different projects. The ideal candidate will be meticulous at the bench, well organized, keep good records and understands the importance of project timelines.

### Job Description

- The successful candidate will spend significant time performing molecular biology bench work including: ATAC-seq, RNA-seq, and RRBS.
- The candidate will record data and assist with analysis.
- The candidate will archive commercial samples, materials and data in a careful and reliable manner.

### Required Skills and Personal Characteristics

- Ability to carefully execute experiments to reliably generate high-quality data according to specified protocols
- An ability to perform and communicate effectively within a group environment
- An ability to learn and implement new molecular biology techniques quickly
- Must possess a strong work ethic and be self-motivated
- Strong interpersonal skills

### Required Experience

- A Bachelor's degree in Biology, Biochemistry, Molecular Biology or other related discipline
- A minimum of one year of relevant laboratory experience with preferred experience in genomics and NGS library preparation
- Proficiency with common software applications (e.g. Excel, Word and PowerPoint)

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](mailto:hr@activemotif.com), referencing "ServicesRA\_2020\_1" in the Subject line.

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