

## RESEARCH ASSOCIATE

Active Motif, Inc. is a privately held biotechnology company focused on supplying innovative kits and reagents for epigenetics and nuclear function research. Active Motif's services group performs specialized epigenetics-based experiments for leading academic groups and pharmaceutical companies and 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 a number of different projects. The ideal candidate will be meticulous at the bench, well organized, keep good records and understand the importance of project timelines.

### Job Description

- The successful candidate will spend significant time performing molecular biology bench work including: chromatin immunoprecipitation (ChIP), qPCR, primer design, and bisulfite sequencing.
- 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

- An ability to carefully execute experiments and 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 self-motivation.

### 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 qPCR and ChIP.
- Proficiency with common software applications (e.g., Excel, Word and PowerPoint).

Interested individuals should submit a cover letter and resume outlining qualifications and experience to [hr@activemotif.com](mailto:hr@activemotif.com).