

**Research Associate, Services (B) – Carlsbad, CA****(Posted: January 4, 2017)**

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 under the supervision of a Research Scientist in the epigenetic services team at our Carlsbad facility. The main job responsibilities will include mastering of Next-Gen sequencing-based techniques such as ChIP-Seq and Reduced Representation Bisulfite Sequencing (RRBS), and generation of data within multiple projects. The ideal candidate will be meticulous at the bench, well organized, keep good records and understand the importance of project time lines.

**Job Description**

The successful candidate will:

- Assist in the development of new techniques
- Perform molecular biology bench work including chromatin immunoprecipitation (ChIP), bisulfite sequencing and high-throughput reporter assays
- Record data and assist with analysis
- 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 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

**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).

Active Motif is an equal opportunity employer and promotes diversity throughout its workforce. Active Motif is headquartered in Carlsbad, California, and has offices in Belgium, Japan and China.

If you are interested in a career with Active Motif, please submit your cover letter and resume outlining qualifications, experience and current salary details to [hr@activemotif.com](mailto:hr@activemotif.com).

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