

Research Associate – Carlsbad, CA

(Posted: June 17, 2016)

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 epigenetics 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 time lines.

The successful candidate will spend significant time performing chromatin preparation using a variety of sample types, chromatin immunoprecipitation (ChIP), qPCR, primer design, bisulfite sequencing and data analysis.

Required Skills

- Ability to follow validated SOPs for sample preparation with close attention to detail
- Ability to carefully execute experiments to reliably generate high-quality data according to specified protocols
- Record data and assist with analysis
- Archive commercial samples, materials and data in a careful and reliable manner
- Contribute ideas & suggestions to improve standard laboratory techniques, improve protocols, processes and equipment
- Perform and communicate effectively within a group environment
- Strong work ethic and 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).

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 located 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.

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