

R&D Associate – Carlsbad, CA

(Posted: Sept. 1, 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 Research and Development group develops assays and kits for use by 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 Research and Development 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 cell culture, protein and DNA extractions and perform protein and DNA-based analysis including chromatin immunoprecipitation and qPCR.

Required Skills

- Ability to carefully execute experiments to reliably generate high-quality data according to specified protocols
- Record data and assist with analysis
- Archive samples, materials and data in a careful and reliable manner
- Contribute ideas and suggestions to improve standard laboratory techniques, 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 a working knowledge of experimental design. Experience in mammalian cell culture, qPCR, epigenetics, and chromatin immunoprecipitation (ChIP) is preferred.
- Demonstrated proficiency in the use of basic laboratory techniques, equipment and materials, and data analysis required
- 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 and Japan.

If you meet the above requirements and are interested in a career with Active Motif, please submit a cover letter and resume outlining qualifications, experience and current salary details to hr@activemotif.com, referencing "R&D Chromatin Associate 2017" in the subject line.

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