

R&D Associate – Carlsbad, CA

(Posted: Sept. 1, 2017)

Active Motif, Inc. is a privately held biotechnology company focused on supplying innovative cell and molecular biology reagents, kits and services for use by leading academic groups and pharmaceutical companies. We are seeking a highly motivated, team-orientated individual to join the R&D team as a Research Associate at our Carlsbad facility.

The Research Associate will work closely with other members of the Research and Development team 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, high throughput cloning, cellular transfections and luciferase assays.

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
- Required laboratory skills include PCR and cloning
- 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 mammalian tissue culture and antibody validation
- 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 Antibody Associate 2017" in the subject line.

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