

R&D Associate II – Antibody Technology (Carlsbad, CA)

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 Research and Development team as a Research Associate at our Carlsbad facility.

The Research Associate will work closely with other members of the R&D 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 timelines. The successful candidate will spend significant time performing cell culture, antibody purification, characterization and validation in a variety of applications (IF, chromatin immunoprecipitation). A minor responsibility of this position is maintenance of laboratory organization and securing competitive quotes for common laboratory supplies.

Required Skills

- Ability to carefully execute, document, and interpret experiments to reliably generate high-quality data according to specified protocols
- Contribute ideas and suggestions to improve standard laboratory techniques, protocols, processes and equipment
- Perform and communicate effectively within a group environment
- Ability to negotiate discounted rates for lab supplies and small equipment
- Required laboratory skills include antibody purification, Westerns, ELISA, IP, and basic DNA techniques including PCR and plasmid purification. Knowledge of IF, FACS or IHC would be an asset.
- Strong work ethic and self-motivated; ability to multitask

Required Experience

- A Bachelor's degree in Biology, Biochemistry, Molecular Biology or other related discipline and three years of relevant laboratory experience, or a Masters' degree and one year of relevant laboratory experience
- Experience in mammalian tissue culture, use of antibodies and biologicals in experiments and PCR
- 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 headquartered in Carlsbad, California, and has additional offices in Belgium, Japan and China.

If you meet the above requirements and are interested in a career with Active Motif, please submit a cover letter and resume outlining your qualifications and experience to: hr@activemotif.com, referencing "R&D Antibody Associate II 2020" in the Subject line.

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