

RESEARCH ASSOCIATE

Active Motif, Inc. is a privately held biotechnology company focused on developing and manufacturing innovative kits and reagents for epigenetics and cell and molecular biology. We are currently looking for a highly motivated, team orientated individual to join our Carlsbad facility as a research associate.

The research associate will work with our R&D and Manufacturing groups to expand our antibody portfolio. The successful individual will be self-motivated, well organized, able to multi-task and be good at time management.

Job Description:

- Plan and design well-controlled experiments.
- Ensure availability of materials and equipment to carry out experiments.
- Make detailed observations, analyze data, and interpret results. Maintain a complete and thorough lab notebook.
- Suggest additional experiments based on research findings.
- Contribute ideas and suggestions to improve standard laboratory techniques, improve protocols, processes and equipment.
- Maintain cell culture facilities and culture cell lines as required for experiments.
- Maintain current lab supplies of common reagents and re-order as needed to maintain workflow.

Requirements:

- Bachelor's degree in Biology, Biochemistry or other related discipline.
- Minimum of one-year laboratory experience, with a working knowledge of experimental design.
- Demonstrated proficiency in the use of basic laboratory techniques, equipment and materials required, tissue culture experience is preferred.
- Experience with Western blots, ELISA, dotblot, immunofluorescence and immunoprecipitation is preferred.
- Experience with hybridoma culturing, antibody production and purification is a plus.

If you are interested, please submit your cover letter and resume outlining your qualifications and experience to hr@activemotif.com. In your cover letter please indicate your experience working with antibodies. We thank all applicants, however, only those selected for interviews will be contacted.