

Research Associate (Carlsbad CA)

Active Motif, Inc. is a privately held biotechnology company that develops, manufactures and supplies innovative cell & molecular biology-based reagents, kits and services that help researchers elucidate the function, regulation and interactions of epigenetic and nuclear processes. We aspire to provide customers with high-quality products, complemented by superior service and support.

We are seeking a highly motivated, team-orientated individual to join the team at our Carlsbad facility. The Research Associate will work in our Manufacturing Department, mostly within the Antibody Production group. The successful candidate will be a highly motivated and detailed oriented researcher with extensive background in cell culture.

Job Description

- Maintenance and expansion of existing and new hybridomas, and manufacturing of monoclonal and recombinant antibodies under manager supervision.
- Maintain cell culture facilities and culture cell lines as required for experiments.
- Follow current protocols and methods and propose new ideas to improve in-house protocols, processes and equipment.
- Make detailed observations, analyze data, and interpret results. Maintain complete and thorough lab notebook.
- Maintain accurate, detailed records, and present updates to managers and project teams.
- Ensure availability of materials and equipment to carry out experiments.
- Ability to multi-task and work on several projects simultaneously.

Requirements

- BS, MS in biological sciences with 1-3 years working experience.
- Impeccable cell culture and sterile technique with demonstrated experience in maintenance and expansion, cryopreservation and cell banking of cancer cell lines a must.
- Experience performing SDS-PAGE required; experience with antibody purification techniques, Western Blotting, ELISA.
- Experience in maintenance and expansion of mouse and rat hybridomas a plus.
- General knowledge of hybridoma methodologies, including immunizations, fusions, selection and screening techniques a plus.
- Experience with medium to large scale production, including bioreactor systems, and recombinant antibody manufacture a plus.
- May be required to work some hours on the weekend for cell maintenance.

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 "Research Associate 2019" in the Subject line. In your cover letter please indicate your experience working with antibodies.

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.

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