

Research Associate (Carlsbad CA)

Active Motif, Inc. is a privately held biotechnology company focused on developing and manufacturing innovative kits and reagents for epigenetics and nuclear function research. We are looking for a highly motivated, team-orientated individual to join us 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 detail-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 team.
- 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 in maintenance and expansion of mouse and rat hybridomas is a plus.
- General knowledge of hybridoma methodologies, including immunizations, fusions, selection and screening techniques is a plus.
- Experience performing SDS-PAGE required; experience with antibody purification techniques, Western Blotting, ELISA.
- Experience with medium- to large scale-production, including bioreactor systems, and recombinant antibody manufacture is a plus.
- May be required to work some hours on the weekend for cell maintenance.

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

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