

Manufacturing 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 currently looking for a highly motivated, team orientated individual to join our manufacturing team to participate in the production and quality control of cell & molecular biology assays at our Carlsbad, CA facility.

Job Description

- Produce buffers, label and aliquot components for assays.
- Run ELISAs, PCR, qPCR, ChIP and Western Blots to validate production quality.
- Maintain cell culture facilities and culture cell lines as required for experiments.
- Follow current protocols and methods as well as 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

- Bachelor's degree in a Biological Science discipline, or the equivalent knowledge and experience.
- A minimum of one-year laboratory experience, with a working knowledge of experimental design. Equivalent experience may be obtained through relevant laboratory class work.
- 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 "Manufacturing Associate, 2019" in the Subject line. In your cover letter, please indicate your experience working with antibodies.

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