

## R&D Research Associate II, Epigenomics / Epitranscriptomics (Carlsbad CA)

Active Motif, Inc. is a privately held biotechnology company focused on developing and supplying innovative products and services for epigenetics and nuclear function research. We are currently looking for a highly motivated, team oriented individual to join the team at our Carlsbad facility.

The Research Associate II will work closely with other members of the Assay team at our Carlsbad facility to develop novel epigenomic / epitranscriptomic workflows for the analysis of complex transcriptional events.

### Job Description

- Develop cutting-edge genome and transcriptome-wide approaches to interrogate DNA and RNA modifications.
- Perform all necessary wet bench experiments for new product development.
- Maintain cultured cell lines (bacteria & eukaryotic) 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 project simultaneously.

### Requirements

- BS or, preferentially, MS in Molecular Biology, Biochemistry, Genomics / transcriptomics, or other related discipline. We will not consider candidates holding a PhD.
- Two - three years laboratory experience.
- Experience with ChIP-Seq; have a detailed understanding of NGS library prep chemistries.
- Demonstrated proficiency in routine molecular biology techniques (e.g. cloning, PCR, qPCR, DNA / RNA extraction).
- Able to work with low-level and nuclease-sensitive analytes such as immunoprecipitated RNA.
- Proficiency in experimental design, data analysis and interpretation.

### Preferences

- Proficient in deep-sequencing techniques such as HiC, GRO-Seq, (h)MedIP-Seq, (h)MerIP-Seq, RIP-Seq, CLIP-Seq.
- Able to run basic bioinformatic analyses.
- Knowledge in DNA/RNA structures such as R-loops, G-quadruplexes, i-Motifs.
- Familiar with lymphoid cell culture (*i.e.* CH12F3-2A or other B cell culture types).
- Use of the Oxford Nanopore for direct DNA and RNA modification analyses.

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.

Interested individuals should submit a cover letter and resume outlining qualifications, experience and availability to [hr@activemotif.com](mailto:hr@activemotif.com), referencing “**Research Associate II**” in the Subject line.

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