

## Associate Bioinformatician (Carlsbad CA)

Active Motif, Inc. is a privately held biotechnology company focused on supplying innovative kits and reagents for epigenetics and nuclear function research. Active Motif's Services group performs specialized epigenetics-based experiments for leading academic groups & pharmaceutical companies and is seeking an Associate Bioinformatician to join our team.

The Associate will work closely with a Senior Bioinformatics Scientist on the Epigenetic Services team to support ChIP-Seq, ATAC-Seq, RNA-Seq and DNA methylation analysis.

### Job Description

- Process raw Illumina sequence data and assign to projects for analysis.
- Run ChIP-Seq, RNA-Seq and ATAC-Seq pipelines, study the underlying code and learn how to customize the analysis.
- Organize and maintain completed project data on fileserver.
- Deliver data and analysis to customers.
- Communicate directly with customers to help them to understand the results.
- Other duties as assigned

### Required Skills and Personal Characteristics

- Highly organized and detail oriented
- Must have experience working from the Linux shell command line
- Knowledge of most commonly used Linux commands, including text editors
- Basic scripting knowledge in Perl, Bash, & R
- Familiarity with Mac OSX and Excel are also important
- Basic understanding of epigenetics and gene regulation

### Required Experience

- A minimum of a Bachelor's degree in biological sciences
- Education and training in the field of bioinformatics
- Some experience with Next-Gen related data analysis

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](mailto:hr@activemotif.com), referencing "Associate Bioinformatician 2019" in the Subject line.

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