

Computational Biologist I (Bioinformatics) (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 and pharmaceutical companies and is seeking a Computational Biologist to join our team.

The Computational Biologist will interact directly with customers and support our epigenetics services and R&D teams at our Carlsbad, CA facility to analyze genome-wide data sets, and report to the Director of Bioinformatics.

Responsibilities

- Provide bioinformatic support to our epigenetics services and R&D team with genome-wide analysis of data sets such as ChIP-Seq, ATAC-Seq, RNA-Seq, HiC, DNA methylation analysis as well as proteomic analysis
- Interact directly with customers to help them understand and interpret results of service projects
- Define and execute analysis pipelines for new services
- Support the improvement and maintenance of current analysis pipelines
- Proactively offer solutions and new computational analysis strategies for novel Epigenomic techniques and data presentation

Required Skills

- Familiarity with Perl, R, and shell scripting
- High attention to detail and scientific & technical problem solving
- Professional written and verbal communication
- Highly organized with maintenance and delivery of large and complex data sets
- Knowledge of Next-Gen sequencing pipelines and Epigenomics

Required Experience

- Ph.D. graduate level training in bioinformatics, computational biology or biology related fields with significant bioinformatics experience
- Experience in Epigenetics and gene regulation
- Expert knowledge of current bioinformatics/genomics/proteomics resources and open source tools

In addition to a highly interactive and stimulating work environment, Active Motif also provides competitive salaries and an excellent benefits package. 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 "Computational Biologist I, Services 2019" in the Subject line.

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