

Technical Support Scientist

Active Motif, Inc. is a privately held biotechnology company focused on supplying innovative kits, reagents and services for epigenetics and gene regulation research. We currently have an opening for a **Technical Support Scientist** to join our team.

Job Description:

We are seeking suitably qualified candidates to provide scientific support and advice to scientists who are using, or are interested in using, Active Motif's high-quality reagents and kits in their research. The selected candidate will provide technical support to scientists throughout the United States and Canada and will collaborate on technical inquiries with our worldwide offices. The role is office-based in Carlsbad, CA, with the potential to attend industry trade shows.

Key Responsibilities:

- Provide technical support to researchers using Active Motif products epigenetics and gene regulation (including ChIP-Seq, DNA methylation analysis and reporter assays).
- Analyze and interpret experimental data, troubleshoot with the customer to resolve the problem.
- Advise customers on the best products and services for their research needs.
- Review and revise new and current product manuals and technical documents
- Writing scientific content for our web pages
- Part of a team involved in new product concept and feasibility.

Required Qualifications:

- Master's or higher degree in Biological Sciences
- 3 or more years laboratory experience in the life science industry
- Enjoy the challenge of addressing diverse scientific problems
- Excellent verbal and written communication skills
- Outstanding organization and time-management skills
- A documented track record of academic and professional success
- Good people skills
- A passion for science!

Desired Qualifications:

- Advanced degree in Epigenetics or Gene Regulation related discipline.
- Good working knowledge of ChIP-Seq, ELISA, Western Blot and Reporter assays
- Previous experience in Technical Support or proven record troubleshooting experiments

Interested individuals should submit a cover letter and resume outlining qualifications and experience to hr@activemotif.com.