

Technical Product Manager (Carlsbad CA)

Active Motif, Inc. is a privately held biotechnology company focused on supplying innovative products and services for epigenetics and nuclear function research. Active Motif is seeking a Technical Product Manager to join our team.

The Technical Product Manager will support our epigenetics services customers and sales team to answer technical questions concerning epigenetic services and develop marketing collateral and campaigns for Epigenetics Services.

Responsibilities

- Interact directly with customers to address technical questions about epigenetic services
- Assist sales team with technical questions about epigenetic services
- Manage, develop, and implement epigenetic services marketing activities
- Maintain sales support documentation
- Perform other duties as required and/or assigned by management

Required Skills

- Professional verbal and written communication skills
- High attention to detail
- Excellent organizational skills
- Team player
- Strong interpersonal skills

Required Experience

- M.S. Biological Sciences or related field
- Marketing experience
- Experience in Epigenetics, Gene Regulation or Molecular Biology
- Working knowledge of Microsoft Office
- Familiarity with Adobe publishing suite and html is preferred
- Customer-facing experience is preferred

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 are interested in a career with Active Motif, please submit your cover letter and resume outlining your qualifications and experience to hr@activemotif.com, referencing “**Technical Product Manager**” in the subject line.

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