

Active Motif, SA - Key Account Manager / The UK (f/m)

Active Motif is the industry leader in developing and delivering innovative tools to enable epigenetics and gene regulation research for the life science, clinical and pharmaceutical and drug discovery communities. We are seeking a highly motivated, independent, and productive team player for the position of Key Account Manager in the UK.

Your tasks

As a member of the Sales team in Active Motif Europe, you will be responsible for the sales of our complete range of our portfolio including Products and Services for epigenetic and gene regulation research throughout the UK. Reporting to the European Sales Manager, your role will involve developing new accounts as well as maintaining existing customers at major academic, biotechnology and pharmaceutical accounts. You will introduce new products and services to our Active Motif customers and support them through in-person meetings, by phone or email. The role will also involve attendance at scientific conferences.

For qualified individuals, the advantages of joining Active Motif include:

- A full and word leading portfolio of assays, genome wide services and validated antibodies for use in ChIP-Seq, backed by thousands of peer-reviewed
- An established customer base, already familiar with Active Motif and our industry-leading product support
- Excellent targeted marketing for rapid expansion into the many labs seeking to begin new epigenetics research projects
- A rich pipeline of innovative technologies, to enable future scientific discovery
- Multidisciplinary Company

Your profile

- You hold a degree Master's or higher degree in Biological Sciences (a PhD and epigenetic background is a plus).
- 3 or more Years laboratory sales experience in the life science industry. (A good knowledge of drug discovery or screening labs is a plus)
- Strong communication and presentation skills
- You enjoy the challenge of addressing diverse scientific problems
- You are an energetic individual who is driven by results
- Outstanding organization and time-management skills
- Regular travel fits your lifestyle (driving license required)
- Based in the UK (London / Cambridge)

Active Motif offers the following rewards:

- Highly competitive salary (based on experience)
- Commission & bonus plans

If you meet these qualifications, and you're excited about an opportunity to join a global leader in epigenetics research tools, forward your CV to: suaud@activemotif.com