

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

---

Active Motif develops & manufactures innovative kits, services and reagent for cell and molecular biology. We seek 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. In addition, you will also support our local distributor with field-based actions and technical support.

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

- A full portfolio of market-leading and innovative epigenetic services, antibodies and research kits, backed by thousands of peer-reviewed publications
- 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

### Your profile

- You hold a degree Master's or higher 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 Cambridge, London or Oxford
- British Citizen

Active Motif offers the following rewards:

- Highly competitive salary (depending 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](mailto:suaud@activemotif.com).