



Fluxion Biosciences is a rapidly-growing cellular assay tool company which is addressing critical needs in the pharmaceutical, biotech, and life science research markets. We are looking for individuals who will thrive in a fast-paced environment and can make meaningful contributions to the sales and commercialization of our line of products for live cell image analysis. Our target assays include automated cell adhesion, microbiological and imaging cytometry. Fluxion is currently looking for a senior level territory manager in the San Francisco Bay area. Responsibilities will entail direct sales and demonstrations of Fluxion products in the Western United States.

Responsibilities:

- Discuss and promote product concepts to prospective customers
- Conduct initial contact to close of sale of Fluxion's products to customer base
- Demonstrate Fluxion products at customer's site
- Provide feedback to product development for product enhancements
- Maintain sales and customer information in corporate database
- Maintain strong relationships with vendors from imaging companies
- Responsible for customer training and installation
- May provide support to other reps outside territory as needed

Requirements

- BS or higher degree in biology-related field
- 5+ yrs selling capital equipment in excess of \$50,000
- A strong rolodex in industry and academia
- Strong communication skills
- Willingness to travel up to 50% overnight within the western United States
- Experience selling into the cell biology arena
- Demonstrates high level of technical and consultative selling skills.
- Preferred candidates will have experience working within a laboratory environment prior to entering sales

Applicants with extensive experience in live cell-based assays will be given first consideration. The successful candidate will become a member of the sales and marketing team and report to the Director of BioFlux Sales. Fluxion offers a competitive compensation package, excellent benefits, and stock options.

Location: South San Francisco, CA

To Apply: Please send Cover Letter and Resume to sales@fluxionbio.com