

Job Title: Research Assistant II (Quality Control)

SignalChem Lifesciences Corporation (SLC) is a clinical stage biotechnology company based in Richmond, British Columbia and is a world recognized leader in kinase research. It consists of a Therapeutics business as well as a Bioreagent business, SignalChem Biotech Inc. (SBI). SBI was established with a focus of producing kinase-related products and other signaling proteins to facilitate the biomedical research and drug discovery efforts of leading research institutions and pharmaceutical companies globally. SignalChem offers signaling protein products encompassing kinases and clinically relevant kinase mutants, phosphatases, epigenetic enzymes, ubiquitin related enzymes, and proteases. SignalChem has established itself as a leader in the development of innovative cell signaling products to meet the needs of the scientist in basic research and drug discovery. The company has utilized its core capabilities of cellular signaling, molecular biology and protein biochemistry to generate around 1500 products as part of its current and continuously expanding product portfolio (currently more than 7000 products). In addition, SignalChem provides research tools and custom services in enzyme activity assay, compound profiling, custom recombinant antibody production and protein production.

In addition, the bioreagent products can also be used to develop diagnostic kits (Biomarkers) to support discovery and development of drug candidates using a personalized medicine approach. This entails identifying those patients with the related kinase defect and applying a specific targeted therapy against the defective kinase which will improve SLC's prospects of successful clinical development and improved patient outcome.

Overview

SignalChem is seeking a well-qualified Research Assistant II to join the Quality Control team. The individual must have exceptional people skills and perform well in a team environment. The position is responsible for planning, conducting, and reporting various experiments in the department of Quality Control.

Duties and Responsibilities

- Sample and test the quality of protein products using current or any associated biochemistry and instrumental techniques by following general guidelines and Standard Operating Procedures (SOPs).
- Make detailed observations, analyze data and interpret results.
- Ensure experimental procedures are carried out carefully and accurately to eliminate errors.
- Prepare, maintain and update databases, technical reports, summaries, and QC protocols.
- Use professional concepts in accordance with QC objectives to solve problems.
- May contribute to the process for creating and executing method development and method validation test protocols for finished products and generating appropriate analytical methods.
- Utilize good laboratory practice including record keeping and lab maintenance.

- Observe and comply with company Health and Safety Policies, Good Manufacturing Practice (GMP) and Good Distribution Practice (GDP).

Requirements

- B.Sc. in Biochemistry or related field with three years of laboratory experience. Comprehensive understanding of quality control.
- Demonstrated track record and skills/experience gained in a similar position(s), at a similar level.
- Highly proficient in protein biochemistry techniques, including SDS-PAGE and the use of key analytical equipment
- Experience with testing, evaluation and release of finished products.
- Self-motivated, with an ability to multi-task and perform under pressure.
- Good computer skills e.g. Microsoft Office (Word, Excel and Outlook). Credible and confident communicator (written and verbal) at all levels.
- Ability to achieve and maintain high standards with meticulous attention to detail with organizational and time management skills.
- Strong analytical and problem solving ability.

Why Work for Us?

SignalChem's employees are passionate, dedicated and extremely motivated to succeed. We are excited by the cutting-edge science and technology, the endless possibilities that we have an opportunity to be a part of. To learn more about SignalChem and our current openings, please visit our website at www.signalchem.com.

We offer challenging career opportunities, competitive benefits and unique work environment.

How to Apply

If you are interested in this challenging opportunity, please send the resumes to humanresources@signalchem.com. Due to the high volume of applicants, only those selected for interviews will be contacted.