

## **Research Associate I**

Luceome Biotechnologies, an emerging biotechnology company, provides unique assay solutions to clients for selectivity profiling of drugs against specific targets. Luceome's assays are based on a technology platform that utilizes luciferase fragment complementation, which allows identification of target and off-target effects of drug candidates with high sensitivity and minimal background interference. For more information, please visit [www.luceome.com](http://www.luceome.com).

## **Position Overview**

Luceome is seeking research oriented, motivated person to contribute to its *KinaseSeeker*<sup>™</sup> program for profiling drug candidates against kinases. The internship candidate will assist in building a large panel of kinases, validating the kinases in split-luciferase based assays, and evaluating test compounds in these assays. Candidates should preferably have research experience in molecular biology techniques, especially cloning, PCR and gel electrophoresis; and biochemical assays. Familiarity with kinases is not required but will be regarded as a plus. Preference will be given to candidates with high analytical ability and excellent communication skills.

## **Qualifications:**

- B.S. or M.S in Biochemistry or Molecular Biology. Candidate should have worked in a laboratory setting for a minimum of 2 years
- Strong track record of scientific achievement
- Hands-on experience in cloning, PCR, gel electrophoresis and biochemical assays
- Willingness to work in a dynamic and changing work environment
- Excellent written and verbal skills

Interested candidates are encouraged to send their CV with cover letter to [info@luceome.com](mailto:info@luceome.com).