



Cell Based Assay Specialist Full-time

Essen Bioscience has an exciting opportunity for a high-calibre, cell-based assay specialist at its European Discovery Services facility on the Biopark, Welwyn Garden City, UK.

The successful applicant will design, assemble, conduct, analyse and interpret cell-based assays to support customer discovery projects and new technology development. This will include new bio-reagent creation (e.g. construction of recombinant cell lines), phenotypic assays using live cell imaging (e.g. angiogenesis, apoptosis) and ion channel assays using planar array automated electrophysiology techniques. Ideally, you will be educated to MSc. level or above in pharmacology, molecular biology, biochemistry or a related discipline, and have >3 years of relevant laboratory experience in a pharmaceutical, biotech or academic environment. Experience in microtitre-plate based assay methods, mammalian cell culture, recombinant techniques, and detailed pharmacological analyses are required. Previous experience with electrophysiology measurements, or in the oncology research field, would be highly valued.

Reporting to the Manager of European Discovery Services, you will be required to analyze, summarize and interpret data for presentation to collaborators and clients. As such you will have strong numeracy, IT, communication and interpersonal skills. You will be a team player, highly motivated, confident and keen to make a difference. Above all, you will be energetic, pro-active and flexible.

In return, we offer a dynamic and stimulating work environment that will challenge and extend your practical and intellectual capabilities. As a growing company, there will be ample opportunity and a strong requirement to learn new scientific and business skills to expand the role over time. A competitive remuneration package including salary, up to 25 days plus bank holidays and free parking will be provided. For further details, or to apply with your CV, contact Jane Moate on 01707 358688 or at moate@essenbio.com (closing deadline 21st August 2011).

Essen Bioscience Ltd.
BioPark, Broadwater Road
Welwyn Garden City
Hertfordshire AL7 3AX, United Kingdom

www.essenbioscience.com