



Ubiquigent Ltd
Sir James Black Centre, College of Life Sciences
University of Dundee, Dow Street
Dundee, Scotland DD1 5EH

July 2015

Job Description

Position: Scientist - Product development and service support

Ubiquigent Ltd

Sir James Black Centre, Dow Street, Dundee, DD1 5EH

Job Summary

The successful applicant will become a critical member of the team in a highly focussed, fast moving, high growth biotechnology product and service company. The role will include supporting and leading the development and delivery of new products and services for customers of Ubiquigent to enable and support their fundamental or drug discovery focused ubiquitin cascade research.

Job Description

Essential skills and laboratory experience:

- Operating and programming liquid handling robotics or other similarly sophisticated automation instrumentation.
- Standard biochemical and cell biology laboratory skills including both *in vitro* and cell culture techniques.
- Experience in the commercial pharmaceutical/biotechnology sectors or in the delivery of services to customers within a non-profit environment.
- Proficiency in the use of Microsoft Excel, Word and Powerpoint.

Desired skills and laboratory experience (2 or more of these skills would be advantageous):

- Knowledge of Microsoft Excel Visual Basic programming.
- Enzymology and/or binding pharmacology.
- Assay development.
- Protein expression and purification; bacteria, baculovirus and/or mammalian cells.
- Product development experience.
- Ubiquitin cascade biology experience.
- Chemical biology research and methods of capturing, purifying and identifying proteins.
- The interaction of the ubiquitin cascade with other cell signalling/post-translational systems; for example phosphorylation.
- The application of advanced analytical techniques – such as MS, ITC, SPR etc – to ubiquitin cascade or other post-translational modification research.



Ubiquigent Ltd
Sir James Black Centre, College of Life Sciences
University of Dundee, Dow Street
Dundee, Scotland DD1 5EH

General requirements:

- A strong drive and ambition to be part of a highly focussed, successful and motivated team in a young high growth biotechnology product and service development company and to have a desire to make a mark and contribute to its continued success and growth.
- Very strong experimental design, data interpretation, problem solving, troubleshooting and communication (both written and oral) skills.
- Willing and able to adapt as may be required when responding to changing circumstances in the scientific field and demands from our customers.
- Desire and ability to learn new techniques and methodologies including the operation of robotics platforms.
- Confident and capable of managing and executing own projects with minimal supervision after development of the ideas and strategy in collaboration with colleagues.
- A strong team player who is motivated both to support and be supported by colleagues.
- A desire to contribute towards the translation of scientific knowledge into products and services to support wider research and understanding of ubiquitin cascade biology and research.
- Motivated and able to contribute original product and service development ideas as well as help refine those of colleagues and collaborators.
- Enthused to be part of a team that collaborates closely with colleagues at our host (University of Dundee) as well as partners from other academic and commercial entities worldwide.
- A strong desire and motivation to provide our customers with the very highest quality products and services and levels of service and support in a timely manner.

Qualifications and Experience

Essential: Relevant HND/HNC degree or BSc or Masters degree and a minimum of three years' experience.

For further details and to apply

Email your CV and a covering letter to careers@ubiquigent.com

Deadline

There is no deadline, however this is a key new role and we're keen to identify the right candidate as soon as possible.



Ubiquigent Ltd
Sir James Black Centre, College of Life Sciences
University of Dundee, Dow Street
Dundee, Scotland DD1 5EH

About Ubiquigent

Ubiquigent was founded in July 2009 and is a high growth biotechnology company whose mission is to be a global leader in developing and supplying the highest quality ubiquitin cascade and related products and services to the life science research and drug discovery markets. The company is venture capital backed and based in the College of Life Sciences at the University of Dundee, Scotland. Products and services are developed and manufactured in-house or developed in collaboration with third parties; both commercial and academic. Ubiquigent has already attracted and delivered projects for major pharmaceutical companies and has an active product and service development programme.

See www.ubiquigent.com.