



Press Release

EsSEN Instruments Announces Name Change and Relocation of Headquarters

Ann Arbor, Michigan – April 12, 2010 – Essen Instruments Inc., a global provider of life science research tools and discovery services, today announced a corporate name change to Essen BioScience Inc. In addition, the Michigan-based company has moved their world headquarters to a newly renovated site at 300 West Morgan Road in Ann Arbor. The new 25,000 sq ft facility provides Essen the space to accommodate the expansion of both their instrumentation and contract service business units.

“Our name change to Essen BioScience better reflects our current global business model” said Kirk Schroeder, President and CEO. “We’ve traditionally been known for innovative life science instrumentation, however our corporate strength is founded in developing novel assays to advance drug discovery. Via our growing service model, we can bring our assays and expertise to more customers”. Commenting further on the corporate move, “We are very proud of our new facility. This investment reflects our commitment to our customers, our employees and to generating jobs in Southeast Michigan.” For further details and contact information visit www.essenbioscience.com.

About Essen BioScience

EsSEN BioScience is a privately held company headquartered in Ann Arbor, Michigan with a wholly-owned subsidiary in Welwyn Garden City, UK. Essen has a long standing tradition of delivering break-through technologies for the life science industry including products such as FLIPR, IonWorks and IncuCyte. To date, their inventions have accounted for nearly 1 billion dollars in worldwide sales. Recently, Essen has expanded their instrumentation-based business model to also include contract services, assembling a team of senior scientists who have successfully managed and driven drug discovery projects at major pharmaceutical companies. The combination of drug discovery and instrumentation experience makes Essen a unique provider of life science research tools and services.

EsSEN BioScience Inc .
300 West Morgan Road
Ann Arbor, Michigan
48108 USA
Tel: 734-769-1600

EsSEN BioScience Ltd.
BioPark, Broadwater Road
Welwyn Garden City, Hertfordshire
AL7 3AX United Kingdom
Tel: 44 (0) 1707 358686