

Media Contacts

D. Doc Chaves, LI-COR Biosciences, dchaves@licor.com, 402-467-0750

Christian Iverson, Quansys Biosciences, civerson@quansysbio.com, 435-752-0531

Quansys Biosciences and LI-COR Biosciences Announce Kits for Infrared Multiplexed ELISA Arrays

*New kits incorporate Quansys' M.E.™ technology
for assays that are central to kinase drug discovery programs*

LOGAN, Utah and LINCOLN, Neb. – Quansys Biosciences™ and LI-COR Biosciences today announced the introduction of four kits for infrared (IR) multiplexed Enzyme Linked Immunosorbant Assay (ELISA) arrays. The kits include Quansys Multiplexed ELISA™ (Quansys M.E.™) Human Cytokine - IR (nine-plex), Human Chemokine - IR (nine-plex), Mouse Cytokine - IR (nine-plex), and Mouse Chemokine - IR (nine-plex). These assays are central to kinase drug discovery programs for indications including oncology, inflammation, metabolic, cardiovascular and neurological diseases.

Quansys Multiplexed ELISA arrays are fully quantitative multiplex ELISA-based tests where up to 16 distinct capture antibodies have been absorbed to each well of a 96-well plate in a defined array. Using less than 30 microliters of sample, up to 84 samples can be assayed for all 16 cytokines in less than two and a half hours, allowing for considerable savings of time, sample and money. The new kits are designed for use on LI-COR Odyssey and Aeries infrared imaging systems.

“We are dedicated to the development of protein arrays to aide researchers and clinicians in better understanding, diagnosing and treating disease for the betterment of people’s lives,” said Matt Groll, vice president of product development for Quansys. “It is our hope that the tools we develop will allow researchers to better serve humanity.”

“These assays add to application flexibility of the Odyssey and Aeries systems and allow the researchers to achieve their research goals on a single platform,” says Neeraj Asundi, LI-COR product marketing manager.

About LI-COR Biosciences LI-COR Biosciences is a leader in the design and manufacture of instrument systems for biotechnology, plant biology and environmental research. The company pioneered the development of infrared fluorescence labeling and detection systems for proteomics, DNA sequencing, microsatellites and AFLP® research. The Odyssey Infrared Imaging System is rapidly becoming the system of choice for quantitative Western Blot analysis. Frost & Sullivan named the Odyssey System the 2006 North American Drug Discovery Technologies Product of the Year.

Founded in 1971, the privately held company is based in Lincoln, Nebraska, with subsidiaries in Germany and the United Kingdom. LI-COR systems are used in more than 100 countries and are supported by a global network of distributors. Further information can be obtained at www.licor.com.

About Quansys Biosciences

Quansys Biosciences, a Spendlove Research Medical Institute (SMRI) company, is dedicated to the development of multiplexed protein arrays to aide researchers and clinicians in better understanding, diagnosing and treating disease. Quansys Biosciences is located in beautiful Cache Valley near Utah State University in Logan, Utah. The SMRI was founded in 1989 and is funded through an endowment established by Dr. Rex S. and Reta Spendlove and family. Dr. Spendlove is the founder of Hyclone Laboratories, a leading producer and distributor of cell culture media throughout the world. Quansys donates a portion of its revenues back to the Institute in an effort to help fund further scientific research that will have a positive impact on human life. Further information can be obtained at www.quansysbio.com