

imaGenes offers first global DNA-Methylation Analysis Service, uses Diagenode MeDIP Kit

Berlin and Liège, February 3, 2009 – imaGenes GmbH, Berlin, adds another unique service – global DNA Methylation Analysis – to its already broad microarray application range. The development of this service offering was greatly assisted by Diagenode’s MagMeDIP Kit, a system using the 5-methyl Cytidine antibody coated on magnetic beads.

DNA methylation is an epigenetic alteration interfering with gene expression. Erroneous DNA methylation may lead to altered gene activity which frequently results in disease. The genome-wide analysis of DNA methylation profiles allows for the identification of novel genetic markers and opens up new avenues for the development of diagnostic and prognostic testing. imaGenes’ comprehensive service offering for the first time enables clinicians as well as fundamental researchers to shed light on the methylation status of patients on a large scale.

“Researchers need to send just 2µg genomic DNA sample to receive extensive DNA methylation profiling data. This removes the requirement for setting-up and optimizing the immunoprecipitation protocol or invest in instrumentation” says Dr. Martina Schad, Associate Director Microarray Services of imaGenes. Raphael Werding, Product Manager of Diagenode concludes “We are very pleased that our High Quality MagMeDIP Kit contributes to the speed and reliability delivered by imaGenes’ MeDIP Service.”

208 words / 1470 characters

About imaGenes :

imaGenes GmbH, Berlin, is a premier provider of genome research services in Europe and was founded in 2007 as a spin-off from RZPD (German Resource Center for Genome Research). imaGenes manages one of the most comprehensive collections of clones of mammalian and vertebrate genes in the world offered via its clone search engine GenomeCube®, linked to international databases. The imaGenes research services feature a broad range of microarray applications on Affymetrix, Agilent, and NimbleGen platforms as well as next generation sequencing services on an Illumina Genome Analyzer System. Researchers in academia and in the pharmaceutical industry are provided with highest quality resources and innovative technology with maximum flexibility and cost effectiveness. imaGenes is fully compliant with ISO 9001:2000 quality standards. A comprehensive overview can be found at www.imagenes-bio.de or www.rzpd.de

Contact:

imaGenes GmbH
Dr. Johannes Maurer – Director Genomic Products & Marketing
Robert-Rössle-Str. 10 / Erwin Negelein Building - 13125 Berlin - Germany
Tel.: +49-30-9489 2440 - eMail: j.maurer@imagenes-bio.de

About Diagenode:

Diagenode sa was founded in 2003 and is the European leading developer and provider of innovative technologies for the field of Epigenetics. The company actively develops a comprehensive portfolio of proprietary products to provide complete solutions from equipments to reagents. Our products are used World Wide by more and more cutting-edge academic and industrial laboratories. These novel products are designed to make assays simpler, faster and more convenient. At the same time we ensure the highest quality results available and the achievement of a much better reproducibility by Automation of these assays. More information about Diagenode or the complete portfolio can be found at www.diagenode.com

Contact:

Diagenode sa
Raphael Werding - Product Manager
Avenue de l'Hôpital, 1 / CHU, Tour GIGA, B34, 3rd floor – Liège - Belgium
Tel.: +32 (0) 4 364 20 56 – eMail: raphael.werding@diagenode.com