



© Alain Janssens

Diagenode is a leading global provider of complete solutions for epigenetics research, biological sample preparation, and diagnostic assays based in Liège, Belgium and Denville, NJ, USA. The company has developed both **(1) shearing solutions for a number of applications, as well as a comprehensive approach to gain new insights into epigenetic studies and (2) innovative products for the infectious diseases market.** The company offers innovative Bioruptor® and Megaruptor® shearing devices, IP-Star® automation instrument, reagent kits, and high quality antibodies to streamline DNA methylation, ChIP, and ChIP-seq workflows. Along with highly trusted and cost-effective In Vitro Diagnostic (IVD) kits for the detection of viruses, bacteria and parasites in a wide number of diseases.

From Diagenode's founding in 2003 in Liège as a local biotechnology startup, the company has expanded rapidly. Diagenode opened its US branch in 2006 and developed a global distribution and partnering network including relationships in Japan and other Asia-Pacific countries.

## 1 EPIGENETICS RESEARCH

Epigenetics plays a revolutionary role in life sciences research. Rapid progress in this field is driving the need for new kinds of research tools and new standards for understanding the epigenetic control mechanisms involved in various biological phenomena such as disease development and drug mechanisms of action. Genome-wide tools which can assess global epigenetic changes at the histone or DNA level (such as Next Generation Sequencing) dramatically advance the understanding of epigenetic mechanisms. Epigenetics research will identify new biomarkers for earlier-stage disease diagnosis and enable development of effective new therapeutics.

Epigenetic assays are complex and necessitate high-quality reagents, robust protocols, and comprehensive data analysis. Diagenode continually innovates solutions and provides the highest quality epigenetics tools available.

## 2 MOLECULAR DIAGNOSTICS

The field of human disease diagnostics is also changing and growing at a rapid rate. Diagenode offers amplification and quantification methods for DNA targets of human infectious disease. We also provide guidance to diagnostics companies developing epigenetics-based diagnostics. Our products are designed to make diagnostic assays simpler, faster, and more convenient, with the highest levels of quality. Our kits are CE-marked and certified ISO 9001 and ISO 13485.

Please contact us for more information [www.diagenode.com](http://www.diagenode.com) [www.diagenodediagnostics.com](http://www.diagenodediagnostics.com)

Europe Diagenode sa / LIEGE SCIENCE PARK / Rue Bois Saint-Jean, 3 / 4102 Seraing (Ougrée) / Belgium / Phone: [+32] 4 364 20 50 / E-mail: [info@diagenode.com](mailto:info@diagenode.com)  
USA Diagenode Inc. / 400 Morris Avenue, Suite 101 / Denville, NJ 07834 / USA / Phone: [+1] 862 209-4680 / Mail: [info.na@diagenode.com](mailto:info.na@diagenode.com)