



Innovating Epigenetics Solutions

## Epigenomics Profiling Services

### Chromatin analysis

- Histone modifications via ChIP-qPCR/ChIP-seq on as low as 20,000 cells/ChIP
- Transcription factor binding via ChIP-qPCR/ChIP-seq
- Chromatin accessibility (bulk ATAC-seq and single cell ATAC-seq)

### RNA sequencing and analysis

- mRNA-seq and gene expression profiling
- Small RNA-seq and miRNA-seq
- Total RNA-seq and whole transcriptome analysis

### DNA Methylation analysis

- Whole genome bisulfite-seq, EPIC array, RRBS, MeDIP/MeDIP-seq
- **Targeted** DNA methylation assays
- **DNA hydroxymethylation** analysis

### Bioinformatic analysis

- Standard analysis including read filtering, mapping and peak calling/methylation calling
- Advanced analysis including differential analysis, annotation, gene ontology, pathway analysis and other customized analysis
- Data mining and machine learning

### Consulting with our scientific experts

- Get expert advice for the best design of your project
- End to end and customized service

	Techniques	Standard analysis	Advanced analysis
Chromatin analysis	<ul style="list-style-type: none"> <li>• ChIP-seq (histones/TFs)</li> <li>• ChIP-qPCR (histones/TFs)</li> </ul>	<ul style="list-style-type: none"> <li>• Read filtering and trimming</li> <li>• Read mapping to the genome</li> <li>• Peak calling</li> </ul>	<ul style="list-style-type: none"> <li>• Identification of differentially bound sites between groups of samples</li> <li>• Genomic annotation</li> <li>• Gene ontology enrichment</li> <li>• Pathway enrichment</li> <li>• Specific genomic region visualization</li> </ul>
	<ul style="list-style-type: none"> <li>• ATAC-seq</li> </ul>	<ul style="list-style-type: none"> <li>• Read filtering and trimming</li> <li>• Read mapping to the genome</li> <li>• Identification of open chromatin regions</li> </ul>	<ul style="list-style-type: none"> <li>• Differentially open chromatin region analysis</li> <li>• Genomic annotation</li> <li>• Nucleosome positioning</li> <li>• Motif recognition</li> <li>• Gene ontology enrichment</li> <li>• Pathway enrichment</li> </ul>
	<ul style="list-style-type: none"> <li>• scATAC-seq</li> </ul>	<ul style="list-style-type: none"> <li>• Read filtering and trimming</li> <li>• Read mapping to the genome</li> <li>• Identification of open chromatin regions</li> <li>• Bead filtration and deconvolution</li> <li>• Cells by peaks count matrix</li> </ul>	<ul style="list-style-type: none"> <li>• Gene ontology enrichment</li> <li>• Pathway enrichment</li> <li>• Cell classification</li> <li>• Differential accessibility</li> </ul>
DNA methylation analysis	<ul style="list-style-type: none"> <li>• RRBS</li> <li>• WGBS</li> </ul>	<ul style="list-style-type: none"> <li>• Read filtering and trimming</li> <li>• Read mapping to the genome</li> <li>• Methylation calling</li> </ul>	<ul style="list-style-type: none"> <li>• CpG and region level differential methylation</li> <li>• Genomic annotation</li> <li>• Gene ontology enrichment</li> <li>• Pathway enrichment</li> </ul>
	<ul style="list-style-type: none"> <li>• MeDIP/hMeDIP-seq</li> </ul>	<ul style="list-style-type: none"> <li>• Read filtering and trimming</li> <li>• Read mapping to the genome</li> <li>• Relative methylation quantification</li> </ul>	<ul style="list-style-type: none"> <li>• Differentially-methylated region identification</li> <li>• Genomic annotation</li> </ul>
	<ul style="list-style-type: none"> <li>• EPIC arrays</li> </ul>	<ul style="list-style-type: none"> <li>• Raw <math>\beta</math>- values</li> <li>• Data normalization</li> <li>• Probe filtering</li> <li>• Background correction</li> </ul>	<ul style="list-style-type: none"> <li>• Gene ontology enrichment</li> <li>• Pathway enrichment</li> </ul>
	<ul style="list-style-type: none"> <li>• Targeted analysis</li> </ul>	<ul style="list-style-type: none"> <li>• Read filtering and trimming</li> <li>• Read mapping to the genome</li> </ul>	<ul style="list-style-type: none"> <li>• Differential methylation</li> </ul>
RNA analysis	<ul style="list-style-type: none"> <li>• RNA-seq</li> </ul>	<ul style="list-style-type: none"> <li>• Read filtering and trimming</li> <li>• Read mapping to the genome or transcriptome</li> <li>• Gene and transcript quantification</li> </ul>	<ul style="list-style-type: none"> <li>• Differentially-expressed gene identification</li> <li>• Annotation</li> <li>• Gene ontology enrichment</li> <li>• Pathway enrichment</li> </ul>
Data mining	<ul style="list-style-type: none"> <li>• Probabilistic modeling</li> <li>• Deep learning approaches</li> <li>• Integrative analyses of different -omics datasets</li> </ul>	<ul style="list-style-type: none"> <li>• Feasibility phase - validation of the data set</li> <li>• Training and testing</li> <li>• Data mining analysis</li> <li>• Interpretation of results</li> </ul>	<ul style="list-style-type: none"> <li>• Reporting</li> <li>• Model validation</li> <li>• Scientific support</li> </ul>

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

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  
USA Diagenode Inc. / 400 Morris Avenue, Suite 101 // Denville, NJ 07834 // USA // Phone: [+1] 862 209-4680 // E-mail: info.na@diagenode.com