

Epigenomics Profiling Services

FROM BIOMARKER DISCOVERY TO BIOMARKER VALIDATION



Consulting with our experts

- Complete end-to-end customized service
- Dedicated scientific team providing high-touch communication
- Expertise with challenging clinical samples



Bioinformatics analysis

- From raw data to differential analysis
- Multi-omics integrative analysis
- Data mining for biomarker discovery



DNA methylation

- Whole genome methylation (WGBS/EM-seq)
- Human Methylome
- RRBS
- Infinium MethylationEPIC v2.0 and Mouse Methylation Array
- Targeted custom methyl-seq



Chromatin

- Histone modifications enrichment or transcription factors binding (ChIP-seq)
- Chromatin accessibility (ATAC-seq)



RNA

- Whole transcriptome (total RNA-seq)
- Messenger RNA (mRNA-seq)