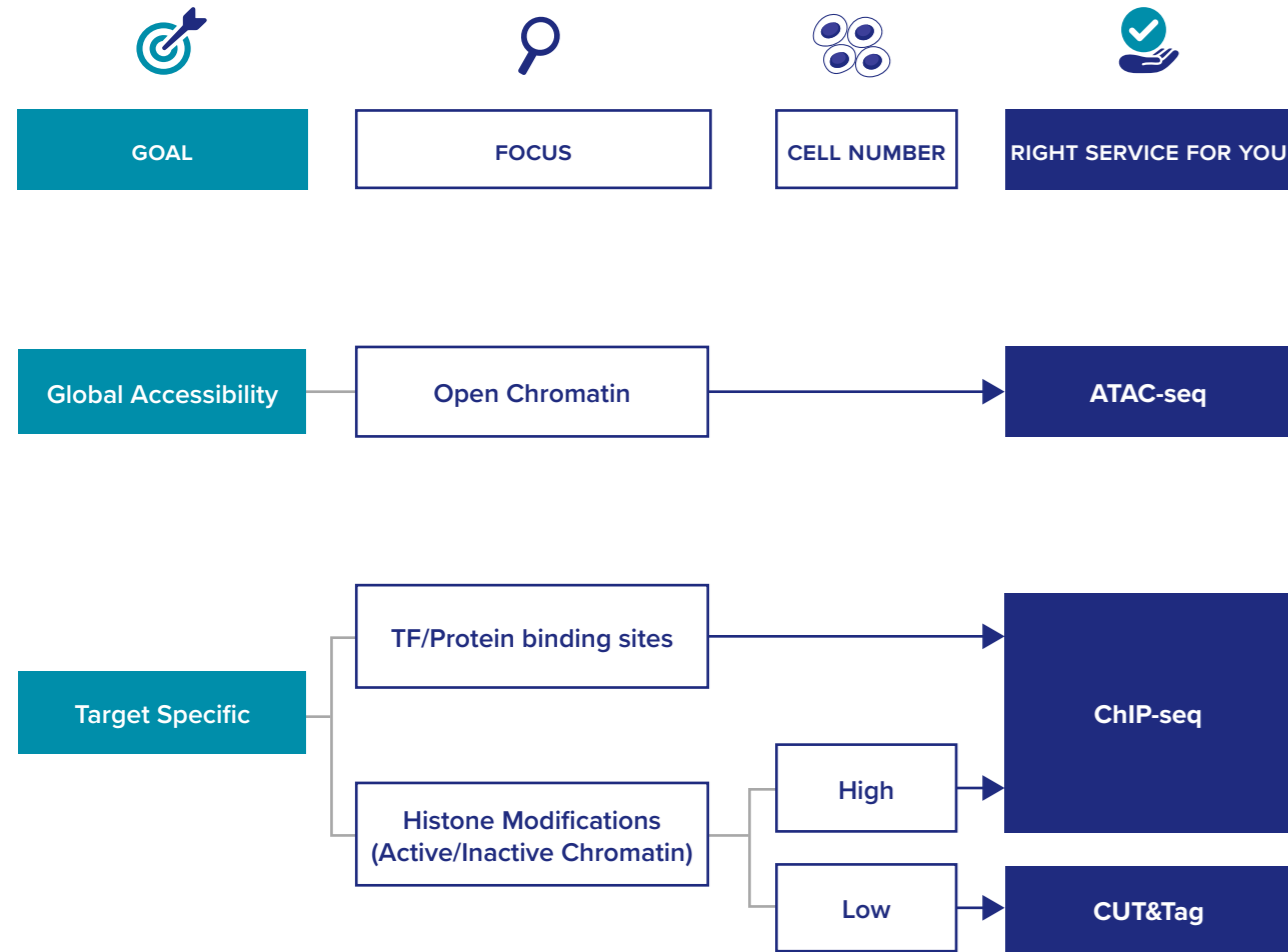
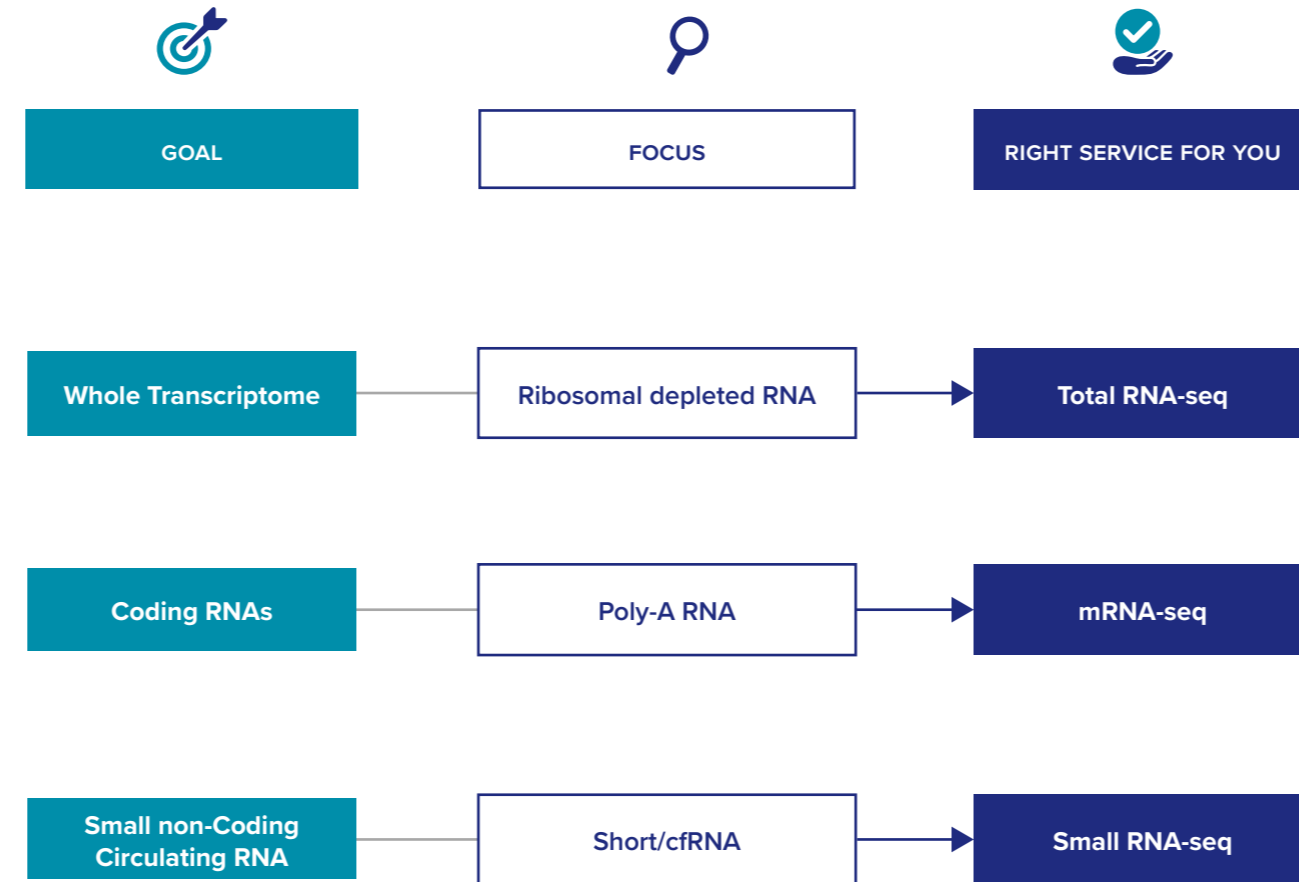


Chromatin



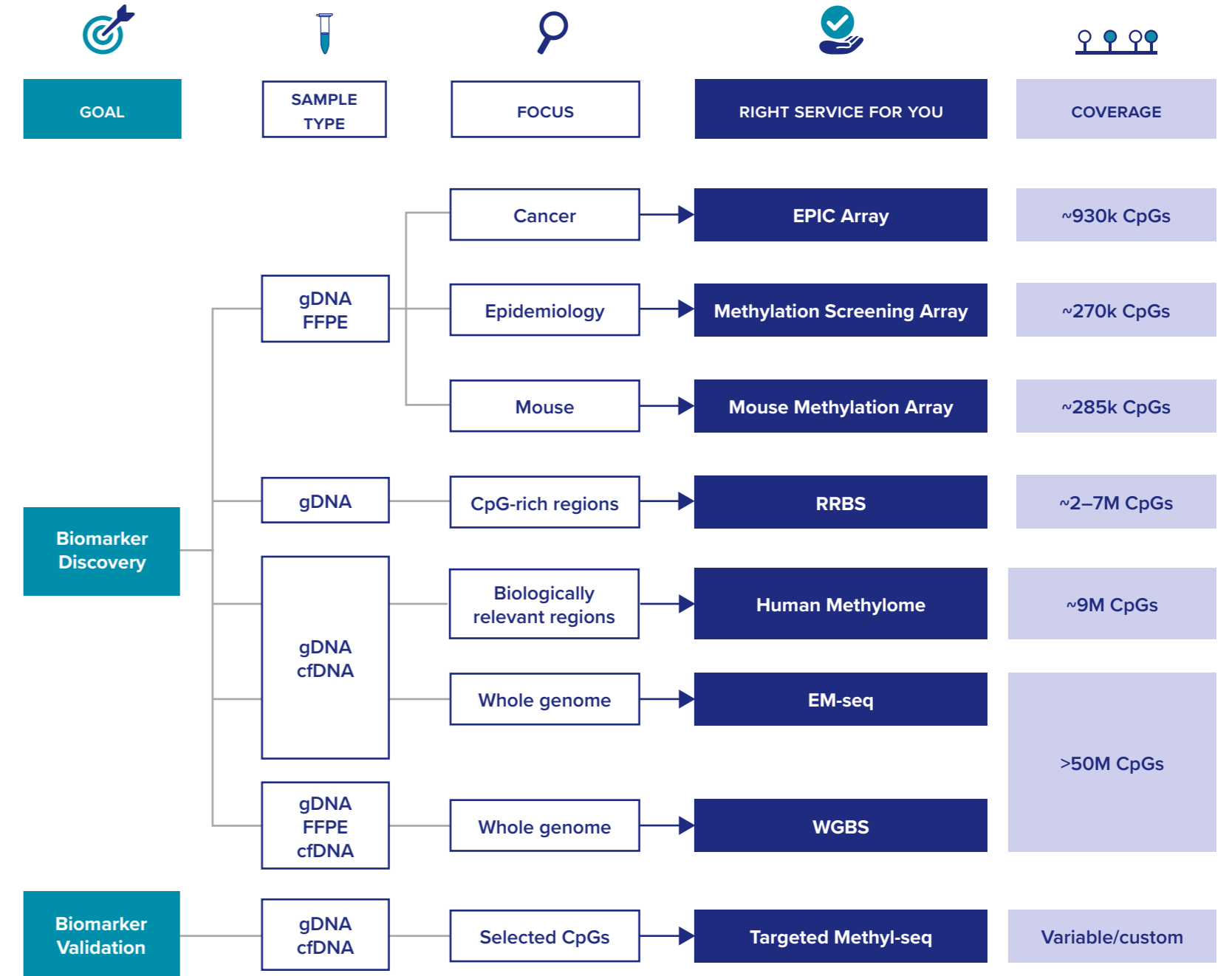
CHROMATIN

RNA



RNA

DNA Methylation



DNA METHYLATION

About Us

Hologic Diagenode stands at the forefront of the epigenomics field, boasting over two decades of expertise. Our advanced services facilitate biomarker discovery and validation by integrating cutting-edge technologies with sophisticated bioinformatics. By delving into the new dimension of epigenomics, we unlock profound insights that transcend genomics, supporting early detection, patient screening, minimal residual disease (MRD) monitoring, and more. We collaborate closely with you to deliver a tailored, end-to-end service, leveraging premier technology to accelerate timelines from R&D to translational research and clinical trials.

Our Service Laboratory Facilities:



USA: San Diego, CA
CLIA environment*

Europe: Liège, Belgium

*Note: Epigenomic testing services are research use only and do not necessarily comply with all CLIA requirements.

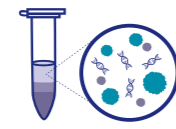
EU // Liège Science Park / Rue Bois Saint-Jean 3, 4102 Seraing / BE / Phone : +32 4 364 20 50 / Fax : +32 4 364 20 51

US // 400 Morris Avenue / Suite 101 Denville / NJ 07834 / USA / Phone : +1 862 209 4680 / Fax : +1 862 209 4681

www.diagenode.com

DNA Methylation in Liquid Biopsies: Overcoming Early Detection Challenges in Cancer

We leveraged our 20 years of epigenomic expertise to help Universal Diagnostics (UDx) develop a customized early-detection workflow for colorectal cancer (CRC), integrating liquid biopsies with DNA methylation technologies for biomarker discovery and validation.



Capturing Biomarkers in Real-Time with Liquid Biopsy

Using cell-free DNA (cfDNA) isolated from plasma, biomarkers were monitored in real-time in the risk population.



DNA Methylation for Cancer Classification

Genome-wide methylation profiles from healthy and cancer patients were created to identify potential biomarkers.



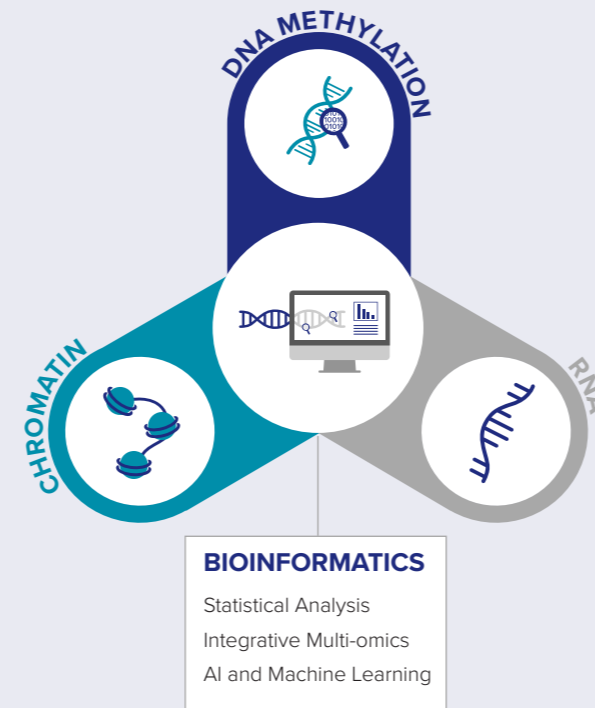
Biomarker Validation

Once a pool of potential biomarkers was available, the best performers were identified.

Learn more



Our Multi-Omics Approach



Our Services by the Numbers



40,000+
samples

including liquid biopsies



250+
projects

Biotech & Pharma



600+
projects

Academia/Government/
Nonprofit Organization (NPO)

AI and Epigenomics: Powering Biomarker Signature Identification

DNA methylation signatures are highly sensitive and specific, making them ideal for early detection, surveillance screening, and monitoring of minimal residual disease (MRD) for cancer and beyond.



DNA Methylation Profiling

Tissue-specific and cancer-specific DNA methylation profiles can be found in previously published and publicly available databases or via genome-wide de novo analysis, such as our WGBS/EM-seq, Human Methylome and EPIC array.



AI and Machine Learning (ML)

Biomarker signatures can be identified bioinformatically using AI and ML analysis approaches.



Biomarker Validation

Biomarkers could then be validated on an independent cohort, publicly available or via new data generated from our custom Targeted Methyl-seq.

Learn more



Why Choose Us?



Profound Biological Insights

Harnessing the power of **epigenomics** combined with advanced **AI** on **clinical samples**, we propel your biomarker discovery to a new dimension.



Accuracy and Reliability

Leveraging our **expertise** in **cutting-edge techniques**, methodologies and platforms, we identify biomarkers with unprecedented precision.



Faster Time to Market

Partner with us to experience a comprehensive journey with **streamlined timelines**, from R&D to clinical trials, accelerating your path to market.



HOLOGIC[®]
Diagenode

Accelerating Biomarker
Discovery with Epigenomics