



**Press Release - May 30, 2013**

## **CDI Laboratories and EpiCypher Form Alliance to Develop Monoclonal Antibodies for Epigenetics Research**

CDI Laboratories and EpiCypher Form Alliance to Develop Monoclonal Antibodies for Epigenetics Research

CDI Laboratories (Mayaguez, Puerto Rico) and EpiCypher, Inc (The Woodlands, TX) have signed an agreement to collaborate on the development of monoclonal antibodies for epigenetics and chromatin biology research. Epigenetics is the study of stable changes in cell function and identity that occur without alterations in DNA sequence.

The project will utilize CDI's Fast-MAb<sup>®</sup> monoclonal antibody development technology and EpiCypher's patented EpiGold<sup>™</sup> Histone Peptide Arrays to develop novel antibodies against histone modifications. Histone modifications are important mediators of epigenetic phenomena, and antibodies to histone modifications are valuable tools in epigenetics research.

"CDI has a terrific setup to develop monoclonal antibodies. Their Fast-MAb<sup>®</sup> system is a perfect fit with our EpiGold<sup>™</sup> Histone Peptide Arrays," stated EpiCypher president Jim Bone.

"EpiCypher's histone peptide array platform allows us to identify antibodies specific for histone modifications quickly and easily," according to CDI vice president and head of business development Scott Paschke. "We are looking forward to a lot of great histone antibodies this year."

About CDI Laboratories: **CDI Laboratories, Inc. (CDI)** is a biotechnology company that delivers products and services founded on whole human proteome microarray technology as well as monoclonal antibody production. CDI's antibodies are tested using its whole human proteome microarray technology, which provides unsurpassed specificity. The company's products and services overcome great hurdles that are currently slowing down advancements in the proteomics field in the areas of research, and protein-based therapeutics and diagnostics.

About EpiCypher: **EpiCypher, Inc.** is a new reagents company founded by renowned chromatin biology researchers to provide novel tools and services for epigenetics and chromatin biology research. EpiCypher's patented EpiGold<sup>™</sup> Histone Peptide Array is the most advanced histone array currently available. EpiCypher also sells nucleosomes, recombinant histone binding proteins, peptides and antibodies, as well as peptide synthesis and peptide array screening services.