



FOR IMMEDIATE RELEASE: 28 January 2014

## **EpiCypher and tebu-bio announce European partnership to distribute industry leading epigenetic products.**

*The pan-European partnership creates a strong combination of the highest quality, innovative products from EpiCypher and the scientific-based relationships of tebu-bio.*

**RESEARCH TRIANGLE PARK, N.C.** – January 28, 2014 – EpiCypher has selected tebu-bio of Le Perray, France to bring European researchers access to EpiCypher’s proprietary products for epigenetics and chromatin biology research. Tebu-bio’s network of local offices in Belgium, France, Germany, Italy, The Netherlands, Denmark, Finland, Norway, Sweden, Portugal, Spain and the U.K. is one of the largest in the European discovery research markets.

Researchers in Europe have a tremendous need for quality products and tebu-bio is a valuable partner to support them. Tebu-bio has supplied European researchers with a diverse portfolio of life science reagents and lab services since 1953. Their commitment to reliable, high quality products is recognized in the market.

“Our partnership with tebu-bio supports our continued commitment to our customers’ success” stated Jim Bone, President of EpiCypher. “During in-person discussions in Europe with tebu-bio, the opportunity to realize our mutual success became clear. As epigenetics gains importance in the global research field, this partnership will increase our presence in the global market to meet demand and further aid the realization of EpiCypher’s business strategy.”

"EpiCypher's portfolio will perfectly complement tebu-bio's offering in the emerging field of epigenetics. On one hand we will explore a new market with EpiCypher's Histone Peptide Arrays: hundreds of highly purified histone derived peptides - unmodified and modified and spotted to the array - allowing in-depth interaction studies”, said Dr. Ali el Baya, tebu-bio Sales Manager for Proteins & Protein based Products & Services. “EpiCypher's native and recombinant nucleosome preparations and epigenetic enzymes will enlarge our offer for screening activities especially in the Drug Discovery environment."

EpiCypher develops and manufactures novel tools for epigenetics and chromatin biology research, based on the expertise of their industry-leading scientific team. Our founders are renowned chromatin researchers, whose knowledge, experience and unique insight into epigenetics benefits our customers in ways that other companies cannot duplicate.

To find out more about EpiCypher’s products available from tebu-bio, visit [tebu-bio's website](#).



### **About EpiCypher**

*Your partner in Epigenetics research.* EpiCypher is a leading epigenetics products and services company that provides research support to leading discovery groups in industry and academia. EpiCypher's patented EpiGold™ 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. For more information about EpiCypher, visit [epicypher.com](http://epicypher.com).

### **About tebu-bio**

tebu-bio is an expert in the supply & distribution of Life Sciences Lab Services & Reagents throughout Europe. Their [own laboratories](#) also provide R&D contract services. Tebu-bio's online catalogue offers Life Science Researchers access to the most innovative range of over 650,000 references. [www.tebu-bio.com](http://www.tebu-bio.com)

### **Contacts:**

Jim Bone  
President  
EpiCypher, Inc.  
855-374-2461  
[jbone@EpiCypher.com](mailto:jbone@EpiCypher.com)

###