



FOR IMMEDIATE RELEASE

EpiCypher Expands Global Distribution Network

Company responds to increased global demand for products and services.

Research Triangle Park, NC - 16 February 2016 - EpiCypher Inc. added five strategic distributors in 2015 responding to increased global demand for the Company's products and services. A total of 16 companies now represent EpiCypher outside of the United States, including nine in Europe.

EpiCypher announced the following distributors after the significant training and support from EpiCypher to enable success in their local markets.

North America:

MJS BioLynx, Inc; Brockville, Ontario, Canada (biolynx.ca)

Asia:

ABclonal Biotechnology Co., Ltd; Wuhan, Hubei, P.R.China (abclonal.com)

Bioinborn Scientific, Inc; Wuhan City, China P.C. (bioinborn.com)

Cistron Biolab, Kattupakkam, Chennai, India (cistronbiolab.com)

KOMA BIOTECH, Seoul, South Korea (komabiotech.com)

"The addition of these highly reputable companies to the EpiCypher distributor portfolio will broaden our sales capacity in established international markets", said Danielle Mahon, Director of Scientific Sales, EpiCypher. "The end-user will benefit from local customer service and expedited shipping of our products."

About EpiCypher - A pioneer in the field of epigenetics and chromatin biology, EpiCypher is a bioscience company developing transformative technologies and delivering superior products to researchers worldwide. EpiCypher's patented EpiTitan™ 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.

Media Contact:

James Bone, Ph.D., President

EpiCypher, Inc.

281-317-7143 (direct)

jbone@epicypher.com

###