

FOR IMMEDIATE RELEASE

EpiCypher Announces Southeast Asian Distribution Channel

Research Triangle Park, NC – 29 March 2021 – Leading epigenetics company, EpiCypher Inc., has added Biomed Global to its world-wide distribution network to strategically address southeast Asia's demand for the Company's products and services. EpiCypher products are available to 27 countries outside of the United States, now including Singapore and Malaysia.

For over 20 years, Biomed Diagnostics Pte Ltd – Biomed Global has been a total solution provider in the biomedical and life sciences industry and is driven by their passion to make a difference to doctors, patients, researchers and academics.

"We're excited to extend distribution rights to Biomed Global. Biomed is a well-established and highly recognized company servicing the life sciences industry in Malaysia and Singapore," said Dr. Jim Bone, President, EpiCypher. "We're confident that their network, local expertise and product knowledge will increase our presence and participation in this important regional market."

Biomed Global (biomed-global.com) represents some of the world's renowned brands known for their products in biomedical sciences, microbiology, molecular biology, genetics testing and imaging systems.

About EpiCypher – EpiCypher was founded in 2012 in response to the growing demand for high-quality reagents and tools to study chromatin regulation and enable epigenetics-focused drug development. The Company develops transformative technologies and offers superior products and discovery services to researchers worldwide. EpiCypher manufactures and sells recombinant "designer" modified nucleosomes (dNucs), including the SNAP-ChIP® product family for quantitative ChIP applications, as well as recombinant histone binding proteins, peptides and antibodies. They also offer a broad range of custom substrate manufacturing and assay development services. Most recently, EpiCypher has been at the leading edge of chromatin mapping technology improvements with the recent launch of highly sensitive epigenomic mapping CUTANA™ assays for ChIC, CUT&RUN, and CUT&Tag.

For more information about EpiCypher, visit **EpiCypher.com**.

To see a complete list of EpiCypher International Distributors, visit **<https://www.epicypher.com/global-distributors/>**.

Media Contact: Leslie Shannon, EpiCypher, Inc. marketing@epicypher.com

###