

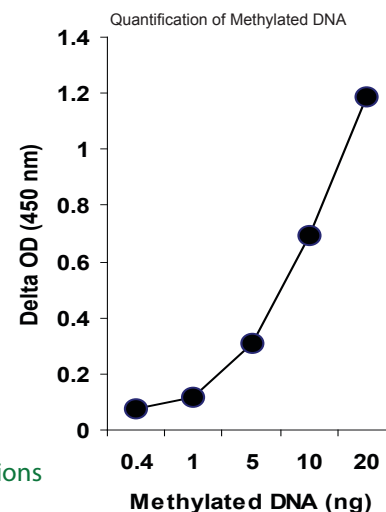
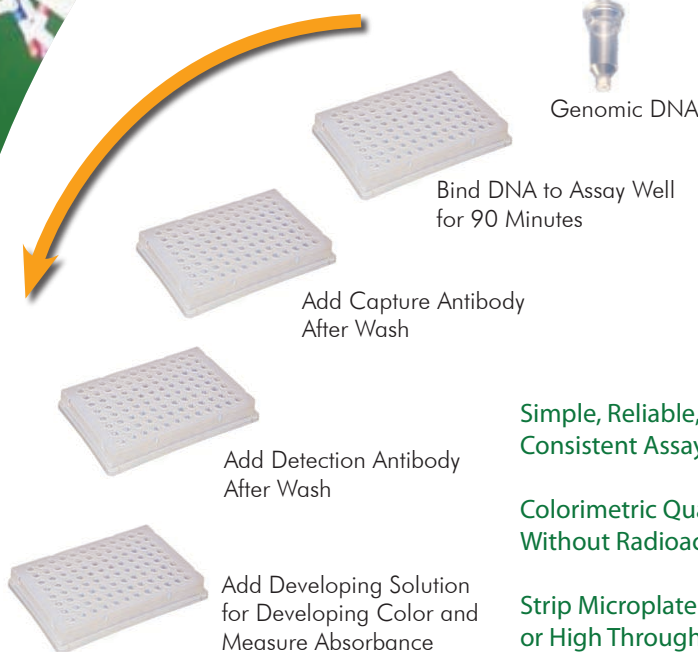
**FINISH IN UNDER
4 HOURS!**

NEW!

Cat No. P-1014B

METHYLAMP™

GLOBAL DNA METHYLATION QUANTIFICATION **ULTRA** KIT



Simple, Reliable, and
Consistent Assay Conditions

Colorimetric Quantification
Without Radioactivity, Extraction, and Chromatography

Strip Microplate Format Makes the Assay Flexible: Manual
or High Throughput Analysis

Rapid Procedure, Which Can Be Finished in
Less Than 4 Hours

High Sensitivity -- Detection Limitation Can
Be as Low as 0.2 ng of Methylated DNA

Universal Positive Control Suitable for Quanti-
fying Methylated DNA From **Any Species**

The Methylamp™ Global DNA Methylation Quantification Ultra Kit is an improved version of the popular Methylamp™ Global DNA Methylation Quantification Kit. As a complete set of essential components, the kit enables the experimenter to quantify global DNA methylation through an ELISA-like reaction using Epigentek's uniquely simplified and streamlined procedure. The entire procedure can be completed in less than 4 hours in either manual or high throughput analysis. In addition to having a shorter procedure, the improved kit also has higher sensitivity than its predecessor. It is suitable for detecting global DNA methylation status using genomic DNA isolated from variety of cultured cells, fresh and frozen tissue, paraffin-embedded tissue, plasma/serum sample, and body fluid sample, etc. The universal positive control included in the kit allows for quantitation of methylated DNA from any species.

PRINCIPLE & PROCEDURE

The Methylamp™ Global DNA Methylation Quantification Ultra Kit contains all reagents required for quantification of global DNA methylation. In this assay, DNA is immobilized to the strip well specifically coated with DNA affinity substance. The methylated fraction of DNA can be recognized by 5-methylcytosine antibody and quantified through an ELISA-like reaction. The amount of methylated DNA is proportional to the OD intensity.

Quantitate Global DNA Methylation Efficiently

ASK YOUR LOCAL DISTRIBUTOR FOR MORE INFORMATION, OR VISIT WWW.EPIGENTEK.COM

Contact Your
Regional Distributor
For Pricing or
Information Today!

Epigentek Group Inc.
United States of America
1 (877) EPIGENTEK
www.Epigentek.com



EPIGENTEK