

April Product of the Month

MethylFlash™ Global DNA Methylation (5-mC) ELISA Easy Kit

*I love how I got results
in just 2 hours!*

DNA Methylation Research
JUST GOT EASIER



- Fast & easy 2-hour ELISA method
- Highly accurate with greater “signal window”
- Detects as low as 0.05% methylated DNA

EPIGENTEK
Complete Solutions for Epigenetics

Take 17% off our MethylFlash Global DNA Methylation
(5-mC) ELISA Easy Kit (Cat. #P-1030)

Use Code: C1704APR

This code is valid April 1 – 30, 2017 and is for US orders only.

Not applicable to distributors or resellers.

www.epigenetek.com · 1-877-374-4368 · sales@epigenetek.com