

January Product of the Month

BisulFlash™ DNA Modification Kit

30-Min Bisulfite Conversion



- Fast 30-min procedure ideal for MSP
- Modified DNA >99.9%
- Input as low as 0.2 ng or 50 cells

EPIGENTEK
Complete Solutions for Epigenetics

Take 17% off our BisulFlash DNA Modification Kit (Cat. #P-1026)*

Use Code: C1701JAN

This code is valid January 1 – 31, 2017 and is for US orders only.
Not applicable to distributors or resellers.

www.epigentek.com · 1877-374-4368 · sales@epigentek.com