

September Product of the Month

ChromaFlash™ High-Sensitivity ChIP Kit

**Insert Chromatin.
Get Results.**



- Easy 5-hour microplate-based method
- Low input of just 2,000 cells
- High specificity with minimal background



Take 17% off our ChromaFlash High-Sensitivity
ChIP Kit (Cat. #P-2027)

Use Code: C1709SEP

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

Not applicable to distributors or resellers.

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