

May Product of the Month

EpiQuik™ Chromatin Accessibility Assay Kit

Open or Closed Detect a Gene's **Chromatin State**



- Finish in only 1 hour & 30 minutes
- Single- or high throughput multi-reaction
- Preserves chromatin structure

EPIGENTEK
Complete Solutions for Epigenetics

Take 17% off our EpiQuik Chromatin Accessibility Assay Kit (Cat. #P-1047)

Use Code: C1705MAY

This code is valid May 1 – 31, 2017 and is for US orders only.

Not applicable to distributors or resellers.

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