

February Product of the Month

EpiQuik™ Histone H3 Modification Multiplex Assay Kit



- 2.5 hour ELISA-like procedure
- 21 well-characterized modifications on 1 plate
- Detects as low as 0.5 ng/well per modification



Take 17% off our EpiQuik Histone H3 Modification Multiplex Assay Kit (Cat. #P-3100)

Use Code: C1702FEB

This code is valid February 1 – 28, 2017 and is for US orders only.

Not applicable to distributors or resellers.

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