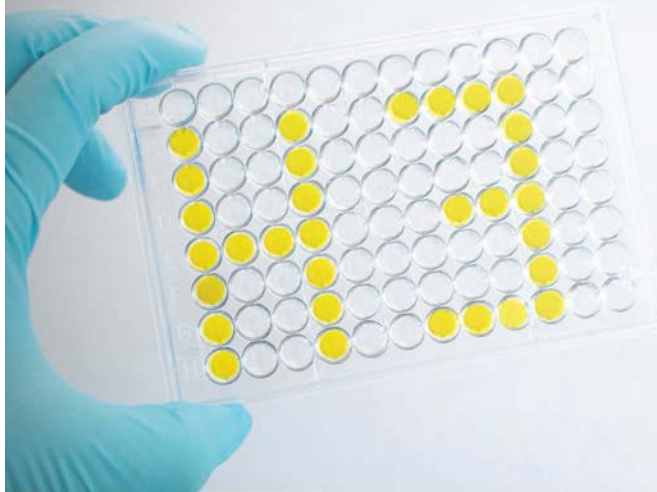


August Product of the Month

Pre-Sure™ Histone H3 Peptide Array ELISA Kit



Quickly Screen for
**Histone H3
Specificity**

- Fast 1 hour & 45 minute procedure
- High quality peptides (over 95% pure)
- Examine 46 different histone modifications



Take 17% off our Pre-Sure Histone H3
Peptide Array ELISA Kit (Cat. #P-3104)

Use Code: C1708AUG

This code is valid August 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