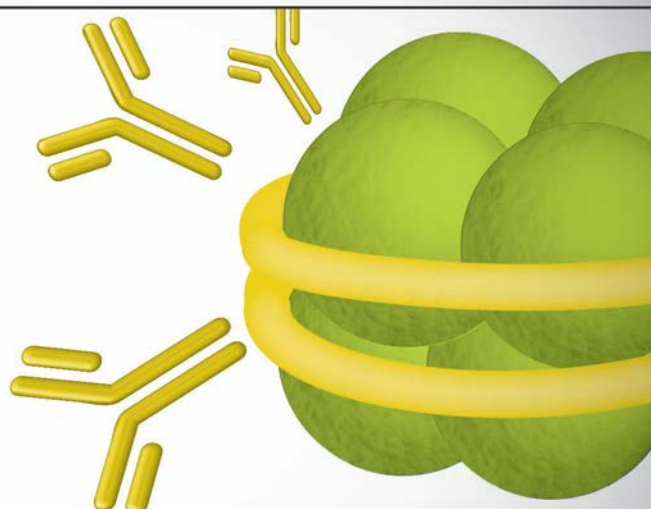


July Product of the Month

Epigenetic Antibodies

**Application
Guaranteed.**



- Specifically for epigenetics & transcription factors
- Risk-free Pioneer Program available for untested species & applications



Take 17% off all our epigenetic antibodies

Use Code: C1707JUL

This code is valid July 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