

## December Product of the Month

### CRISPR/Cas9 Monoclonal Antibody



for precise gene  
editing studies

- Detect CRISPR/Cas9 expression levels to confirm successful transfection
- Detect genome-wide effects of CRISPR/Cas9 by Cas9 ChIP



Take 17% off our CRISPR/Cas9 Monoclonal  
Antibody [7A9] (Cat. #A-9000)

**Use Code: C1712DEC**

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

Not applicable to distributors or resellers.

[www.epigentek.com](http://www.epigentek.com) · 1-877-374-4368 · [sales@epigentek.com](mailto:sales@epigentek.com)