

March Product of the Month

EpiQuik™ m6A RNA Methylation Quantification Kit

Reveal the Mysteries of RNA Methylation



- Simple, 4-hour ELISA-like procedure
- Detects as low as 10 pg of m6A
- Quantify m6A from mRNA & ncRNA



Take 17% off our EpiQuik m6A RNA Methylation
Quantification Kit (Cat. #P-9005)

Use Code: C1703MAR

This code is valid March 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