

## June Product of the Month

### EpiQuik™ Circulating Cell-Free DNA (cfDNA) Isolation Easy Kit

Magnetic Bead-Based Purification of  
**Small Fragment  
Circulating cfDNA**  
from Plasma or Serum



- Fast, straightforward 1-hour procedure
- DNA suitable for qPCR & NGS DNA library prep
- High recovery >80% input DNA

**EPIGENTEK**  
Complete Solutions for Epigenetics

Take 17% off our EpiQuik Circulating Cell-Free DNA (cfDNA)  
Isolation Easy Kit (Cat. #P-1065)

**Use Code: C1706JUN**

This code is valid June 1 – 30, 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)