

November Product of the Month

EpiQuik™ 8-OHdG DNA Damage Quantification Direct Kit



- Less than 4 hour ELISA-like procedure
- Detection limit as low as 2 pg of 8-OHdG
- No cross-reactivity to 8-OHdG analogues



Take 17% off our EpiQuik 8-OHdG DNA Damage
Quantification Direct Kit (Cat. #P-6003)

Use Code: C1711NOV

This code is valid November 1 – 30, 2017 and is for US orders only.

Not applicable to distributors or resellers.

www.epigentek.com · 1-877-374-4368 · sales@epigentek.com