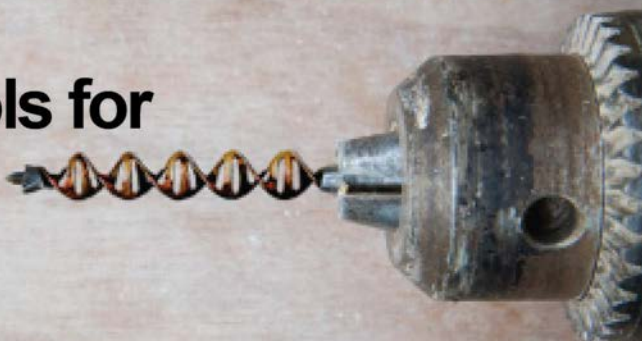


October Product of the Month

Epigenase™ 5mC-Hydroxylase TET Activity/Inhibition Assay Kit

We Have the Tools for
TET Studies



- Fast, easy-to-follow 5-hour procedure
- Directly measures TET hydroxylase activity
- Extremely low background signals



Take 17% off our Epigenase 5mC-Hydroxylase
TET Activity/Inhibition Assay Kit (Cat. #P-3086)

Use Code: C1710OCT

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

Not applicable to distributors or resellers.

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