

## Epigenetic Detection of Activity of HMT in Cancer Cells

### (Measurement of Methylation Modification in Cells)

**Cat #: HMT-Western**

**Cells: Human Prostate Cancer Cells: DU145, PC-3**

#### **Description:**

Human Prostate cancer cell lines, DU145 and PC-3, were treated with 10 $\mu$ M UNC0638 (U), 5 $\mu$ M SAHA (S), 1 $\mu$ M TSA (T) or DMSO (D) control for 2 days in the presence of 10% fetal bovine serum. Whole cell lysates of the treated cells were subject to Western blot analyses with anti-dimethyl-Histone H3K9 antibody or control antibody anti- $\alpha$ -Tubulin.

