

Epigenetic Detection of Activity of HDAC in Cancer Cells

(Measurement of Acetylation Modification in Cells)

Cat #: HDAC-Western

Cells: Human Prostate Cancer Cells: DU145, PC-3; Human Fibroblast, MRC5

Description:

Human fibroblast cell line, MRC5, and human prostate cancer cell lines, DU145 and PC-3, were treated with 10 μ M Tubastatin A (Tu), 5 μ M SAHA (S), 1 μ M TSA (Ts), or DMSO (D) as control in the presence of 10% FBS for 22 hrs. Whole cell lysates of the treated cells were subject to Western blot analyses with anti-Acetylated α -Tubulin, anti- α -Tubulin, anti-acetylated Histone H3 or anti-Histone H3 antibodies.

