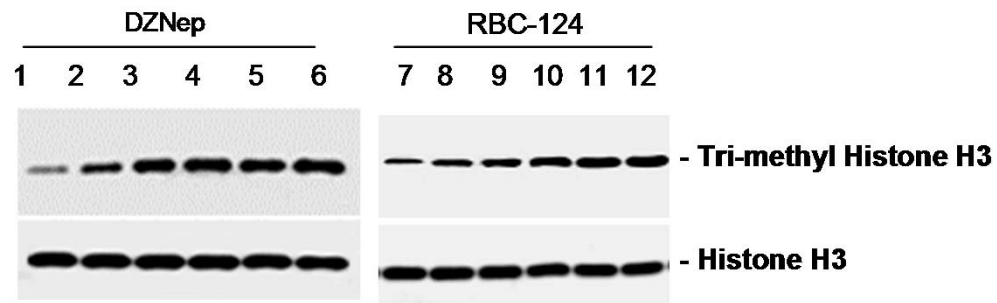


Measurement of Histone H3 Methylation in Cells by Western Blot

**Tri-methylated or Total Histone H3 in SUDHL-6 cells
treatment with compound and measure by Western Blot**



Sample #	DZNep (uM)	Sample #	RBC-124 (uM)
1	10.000	7	100.00
2	2.000	8	20.00
3	0.400	9	4.00
4	0.080	10	0.80
5	0.016	11	0.16
6	DMSO ctrl	12	DMSO ctrl