

Measurement of H3K27ME3 Level in cells using AlphaLISA Format

Su-DHL-6 Cells: Human B cell lymphoma cell line from ATCC.

Assay format: Cell-based AlphaLISA assay.

Assay Protocol: 5000 SU-DHL-6 cells were seeded to assay plates. Cells were incubated with compound for 3 days at 37°C in 5% CO₂. Evaluation of the cellular inhibitory activity of the reference inhibitor generated an average IC₅₀ value of 13.0 ± 1.3 μM and 1.4 ± 0.4 μM for RBC124 (■) and DZNep (▼), respectively. The IC₅₀ (EC₅₀) values were calculated and the IC₅₀ curves were plotted using the GraphPad Prism 4 program based on a sigmoidal dose-response equation.

