

NanoBRET™ Target Engagement Assay for EPHA5 Kinase

Cell line: HEK293 human epithelial cell line from ATCC.

Assay format: 384-well plate format using Promega's NanoBRET™ TE Intracellular Kinase Assay.

Assay protocol: HEK293 cells transiently expressing NanoLuc®-EPHA5 Fusion Vector were seeded into the wells of 384-well plates. The cells were pre-treated with the NanoBRET™ Tracer K-4 and then treated with reference compound Dasatinib for 1 hour. The BRET signal was measured on an Envision 2104 Multilabel Reader. IC₅₀ value was calculated and IC₅₀ curve was plotted using the GraphPad Prism 4 program based on a sigmoidal dose response equation.

Reference compound: Dasatinib (IC₅₀ value): 1.303e-009 M

