

NanoBRET™ Target Engagement Assay for MAP4K2 Kinase

Cell line: HEK293 human epithelial cell line from ATCC.

Assay format: 384-well plate format.

Assay protocol: HEK293 cells transiently expressing MAP4K2-NanoLuc fusion vector were seeded into the wells of 384-well plates. The cells were pre-treated with the tracer 05 and then treated with reference compound CTx-0294885 for 1 hour. NanoBRET™ signal was measured on an Envision 2104 Multilabel Reader. IC50 (EC50) value was calculated and IC50 curve was plotted using the GraphPad Prism 4 program based on a sigmoidal dose-response equation.

Reference compound: CTx-0294885 (IC50 value): 8.446e-007 M

