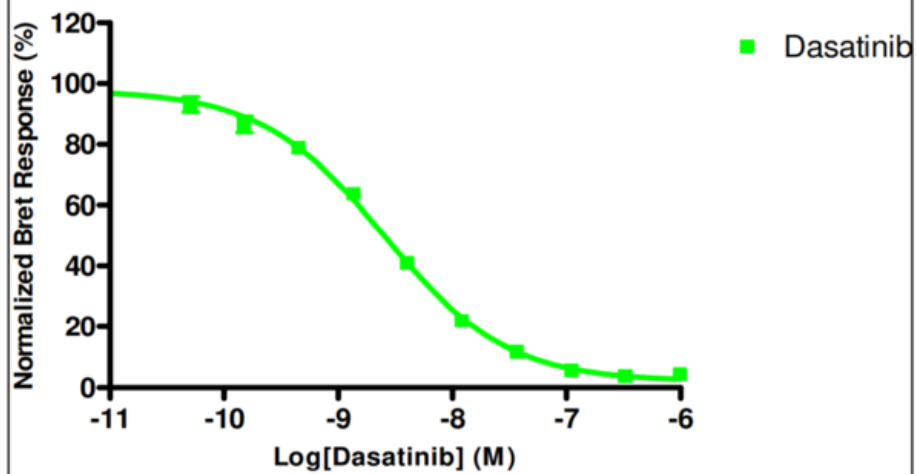


EPHA2

NanoBRET Target Engagement Intracellular Kinase Assay, Competitive binding assay format

Target	EPHA2				
HGNC Symbol	EPHA2				
Synonyms	ECK				
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-EPHA2 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.				
Readout	Binding of compound to target kinase				
Reference Compound IC50s	<table border="1"> <thead> <tr> <th>Compounds</th> <th>IC50 (M)</th> </tr> </thead> <tbody> <tr> <td>Dasatinib</td> <td>2.528e-009</td> </tr> </tbody> </table>	Compounds	IC50 (M)	Dasatinib	2.528e-009
Compounds	IC50 (M)				
Dasatinib	2.528e-009				
Screening Location	Malvern, USA				

Reference Compound IC50 for EPHA2



HILLSLOPE	-0.8109
EC50	2.528e-009