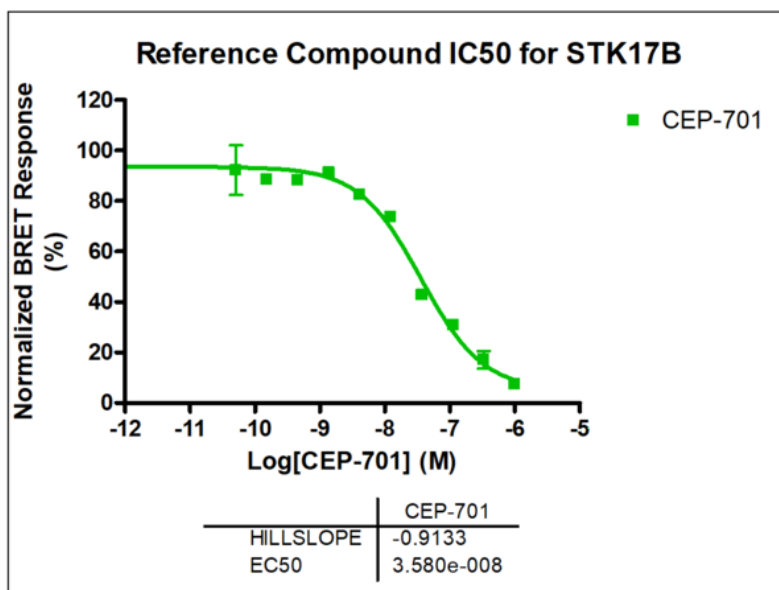


STK17B NanoBRET Kinase Assay

STK17B [NanoBRET Target Engagement Intracellular Kinase Assay](#) for screening of compounds binding to the target.

Target	STK17B				
HGNC Symbol	STK17B				
Synonyms	DRAK2				
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- STK17B Fusion Vector were seeded into the wells of 384-well plates. The cells were pre-treated with the NanoBRET Tracer K-9 and then treated with reference compound CEP-701 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>CEP-701</td> <td>3.580e-008</td> </tr> </tbody> </table>	Compounds	IC50 (M)	CEP-701	3.580e-008
Compounds	IC50 (M)				
CEP-701	3.580e-008				
Screening Location	Malvern, USA				



Let's Discover Together.

