

# SIK2

## NanoBRET Target Engagement Intracellular Kinase Assay, Competitive binding assay format

<b>Target</b>	SIK2				
<b>HGNC Symbol</b>	SIK2				
<b>Synonyms</b>	SIK2				
<b>Cell Line</b>	HEK293 human epithelial cell line from ATCC				
<b>Assay Format</b>	384-well plate format using Promega's NanoBRET TE Intracellular Kinase Assay				
<b>Assay Protocol</b>	HEK293 cells transiently expressing NanoLuc-SIK2 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.				
<b>Readout</b>	Binding of compound to target kinase				
<b>Reference Compound IC50s</b>	<table border="1"> <thead> <tr> <th>Compounds</th> <th>IC50 (M)</th> </tr> </thead> <tbody> <tr> <td>Dasatinib</td> <td>1.447e-008</td> </tr> </tbody> </table>	Compounds	IC50 (M)	Dasatinib	1.447e-008
Compounds	IC50 (M)				
Dasatinib	1.447e-008				
<b>Screening Location</b>	Malvern, USA				

### Reference Compound IC50 for SIK2

