

# MAP4K5 NanoBRET Kinase Assay

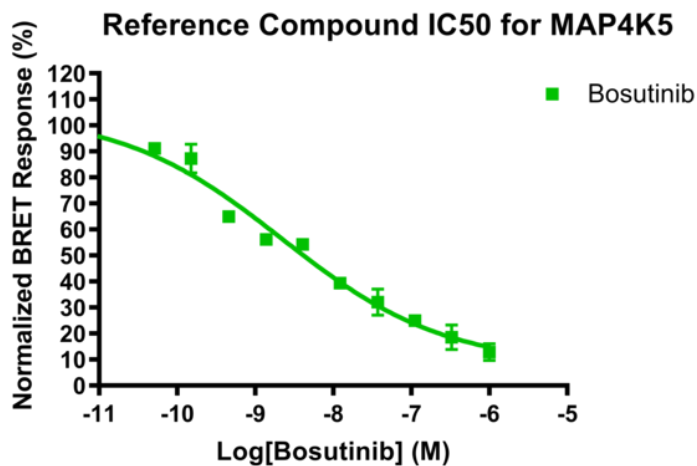
MAP4K5 [NanoBRET Target Engagement Intracellular Kinase Assay](#) for screening of compounds binding to the target.

## Target Overview

Target	MAP4K5
HGNC Symbol	MAP4K5
Synonyms	GCKR, KHS, KHS1, MAPKKKK5
Cell Line	HEK293 human epithelial cell line from ATCC

## Assay Properties

Assay Format	384-well plate format using Promega's NanoBRET TE Intracellular Kinase Assay	
Assay Protocol	HEK293 cells transiently expressing NanoLuc- MAP4K5 Fusion Vector were seeded into the wells of 384-well plates. The cells were pre-treated with the NanoBRET Tracer K-10 and then treated with reference compound Bosutinib 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	<b>Compounds</b>	<b>IC50 (M)</b>
	Bosutinib	2.218e-009
Screening Location	Malvern, USA	



	Bosutinib (M)
HillSlope	-0.4168
EC50	2.218e-009

Let's Discover Together.

© Oct 2021 Reaction Biology. All rights reserved.

[https://www.reactionbiology.com/datasheet/map4k5\\_nano\\_malvern](https://www.reactionbiology.com/datasheet/map4k5_nano_malvern)

