

# KRAS (G12V) NanoBRET Ras Assay

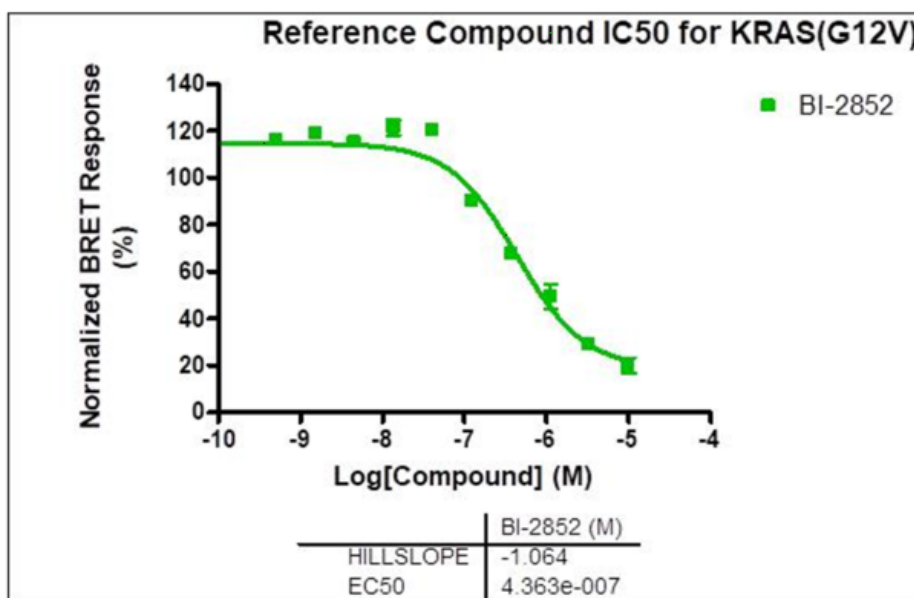
KRAS (G12V) [NanoBRET Target Engagement Intracellular Ras Assay](#) for screening of compounds binding to the target.

## Target Overview

Target	KRAS (G12V)
HGNC Symbol	KRAS
Synonyms	KRAS2B, KRAS2A, KRAS4A, KRAS4B, RASK2
Cell Line	HEK293 human epithelial cell line from ATCC

## Assay Properties

Assay Format	384-well plate format using Promega's NanoBRET TE Intracellular Ras Assay	
Assay Protocol	HEK293 cells transiently expressing LgBiT®-KRAS (G12V) and SmBiT®-KRAS (G12V) Fusion Vectors were seeded into the wells of 384-well plates. The cells were pre-treated with the NanoBRET Tracer K-1 and then treated with reference compound BI-2852 for 2 hours. 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 Ras	
Reference Compound IC50s	<b>Compounds</b>	<b>IC50 (M)</b>
	BI-2852	4.363e-007
Screening Location	Malvern, USA	



Let's Discover Together.

