

# ACVR1 (Q207D) NanoBRET Kinase Assay

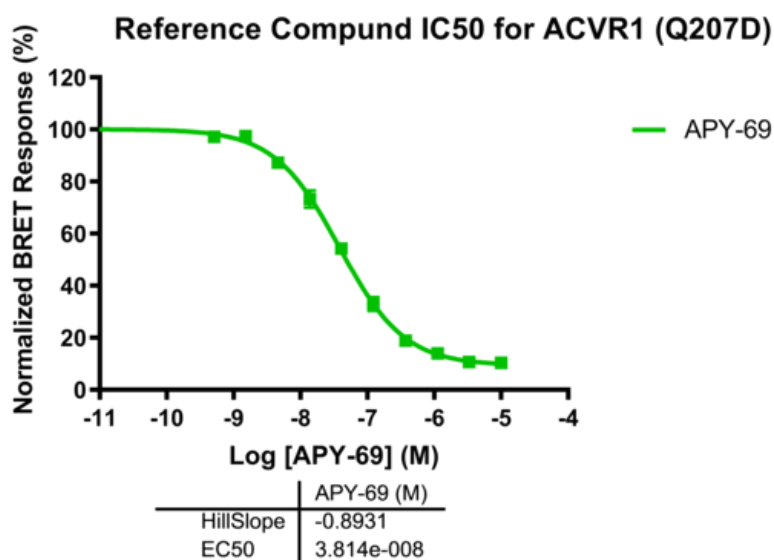
ACVR1 (Q207D) [NanoBRET Target Engagement Intracellular Kinase Assay](#) for screening of compounds binding to the target.

## Target Overview

Target	ACVR1 (Q207D)
HGNC Symbol	ACVR1 (Q207D)
Synonyms	ACTRI, ACVR1A, ACVRLK2, ALK2, FOP, SKR1
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- ACVR1 (Q207D) Fusion Vector were seeded into the wells of 384-well plates. The cells were pre-treated with the NanoBRET Tracer K-11 and then treated with reference compound APY-69 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	<b>Compounds</b>	<b>IC50 (M)</b>
Compound IC50s	APY-69	3.814e-008
Screening Location	Malvern, USA	



Let's Discover Together.

