

# BRD4-2 NanoBRET TE Intracellular Assay

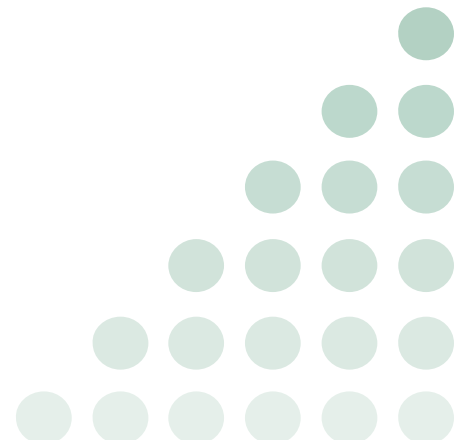
BRD4-2 [NanoBRET Target Engagement Intracellular Assay](#) for screening of compounds binding to the target, competitive binding assay format

## Target Overview

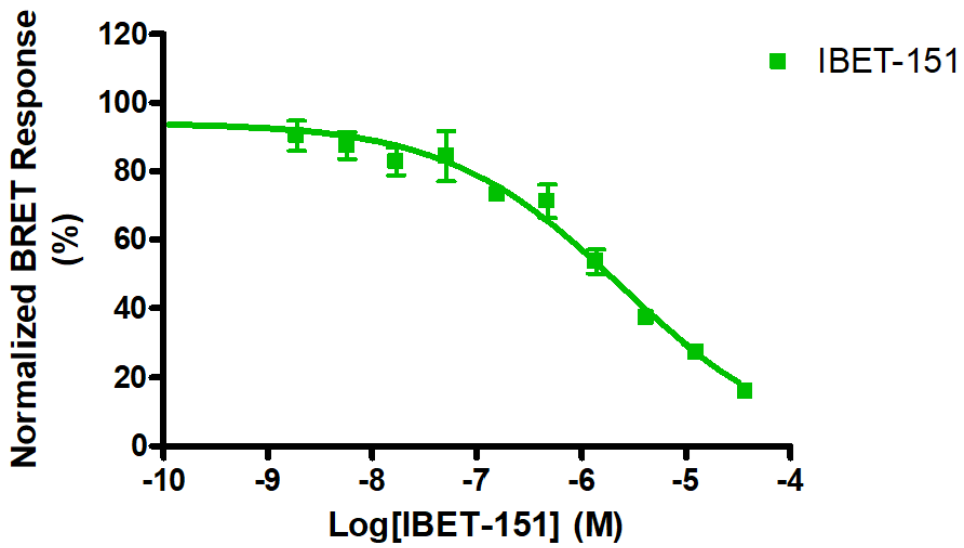
<b>Target</b>	BRD4-2
<b>HGNC Symbol</b>	BRD4 (BD2)
<b>Synonyms</b>	HUNK1; MCAP; CAP; HUNK1
<b>Cell Line</b>	HEK293 human epithelial cell line from ATCC

## Assay Properties

<b>Assay Format</b>	384-well plate format using Promega's NanoBRET Intracellular Target Engagement Assay
<b>Assay Protocol</b>	HEK293 cells transiently expressing NanoLuc-BRD4 (BD2) Fusion Vector were seeded into the wells of 384-well plates. The cells were pre-treated with the NanoBRET BRD Tracer and then treated with reference compound IBET-151 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 program based on a sigmoidal dose response equation.
<b>Readout</b>	Binding of compound to target
<b>Reference Compound IC50s</b>	<b>Compounds IC50 (M)</b> IBET-151 2.257e-006
<b>Screening Location</b>	Malvern, PA, USA
<b>Further information</b>	More information can be found on our website <a href="#">Cell-based Epigenetic Assays</a>



Reference compound IC50 for BRD4 (BD2)



	IBET-151 (M)
HILLSLOPE	-0.5272
EC50	2.257e-006

