

MAP3K3 NanoBRET Kinase Assay

MAP3K3 [NanoBRET Target Engagement Intracellular Kinase Assay](#), competitive binding assay format for screening of compounds binding to the target.

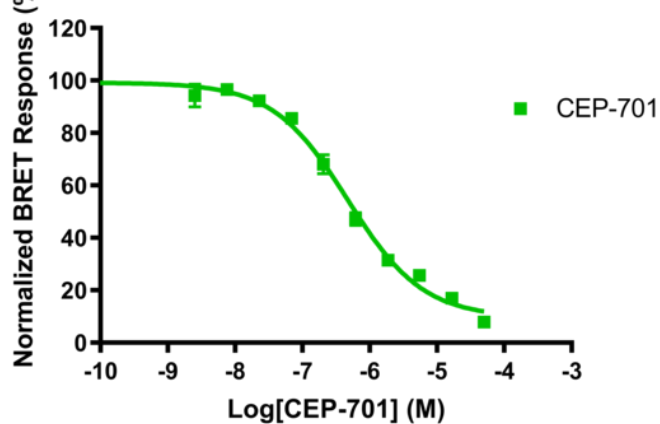
Target Overview

Target	MAP3K3
HGNC Symbol	MAP3K3
Synonyms	MAPKKK3, MEK kinase 3, MEKK3
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-MAP3K3 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 CEP-701 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	Compounds IC50 (M) CEP-701 4.7330e-007
Screening Location	Malvern, USA

Reference Compound IC50 for MAP3K3



Let's Discover Together.

© Dec 2022 **Reaction Biology**. All rights reserved.

https://www.reactionbiology.com/datasheet/map3k3_nano_malvern

