

# PIK3CA (E545A)/PIK3R1 NanoBRET Kinase Assay

PIK3CA (E545A)/PIK3R1 [NanoBRET Target Engagement Intracellular Kinase Assay](#) for screening of compounds binding to the target.

Target Overview					
Target	PIK3CA (E545A)/PIK3R1				
HGNC Symbol	PIK3CA (E545A)-PIK3R1				
Synonym	MGC142161, MGC142163, p110-alpha, PI3K, PI3K-alpha				
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 co-expressing NanoLuc-PIK3CA (E545A) Fusion Vector and PIK3R1 Expression Vector were seeded into the wells of 384-well plates. The cells were pre-treated with the NanoBRET Tracer K-3 and then treated with reference compound PI-103 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.				
Readout	Binding of compound to target kinase				
Reference Compound IC50s	<table border="1"> <thead> <tr> <th>Compound</th> <th>IC50 (M)</th> </tr> </thead> <tbody> <tr> <td>PI-103</td> <td>3.932E-008</td> </tr> </tbody> </table>	Compound	IC50 (M)	PI-103	3.932E-008
Compound	IC50 (M)				
PI-103	3.932E-008				
Screening Location	Malvern, PA, USA				
Further Information	More information can be found on our website <a href="#">NanoBRET Target Engagement Intracellular Kinase Assay Services</a> .				
Reference compound IC50 for PIK3CA (E545A)/PIK3R1					

Let's Discover Together.

© May 2024. Reaction Biology. All rights reserved.

