

ULK3 NanoBRET Kinase Assay

ULK3 [NanoBRET Target Engagement Intracellular Kinase Assay](#) for screening of compounds binding to the target.

Target Overview					
Target	ULK3				
HGNC Symbol	ULK3				
Synonym	FLJ90566, DKFZP434C131				
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-ULK3 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 Staurosporine 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>Staurosporine</td> <td>3.048E-007</td> </tr> </tbody> </table>	Compound	IC50 (M)	Staurosporine	3.048E-007
Compound	IC50 (M)				
Staurosporine	3.048E-007				
Screening Location	Malvern, PA, USA				
Further Information	More information can be found on our website NanoBRET Target Engagement Intracellular Kinase Assay Services .				
Reference compound IC50 for ULK3					

Let's Discover Together.

© May 2024. Reaction Biology. All rights reserved.

