

IGF1R Kinase Assay Service

Activity-based IGF1R kinase assay for compound screening and profiling via [radiometric HotSpot™ kinase assay](#).

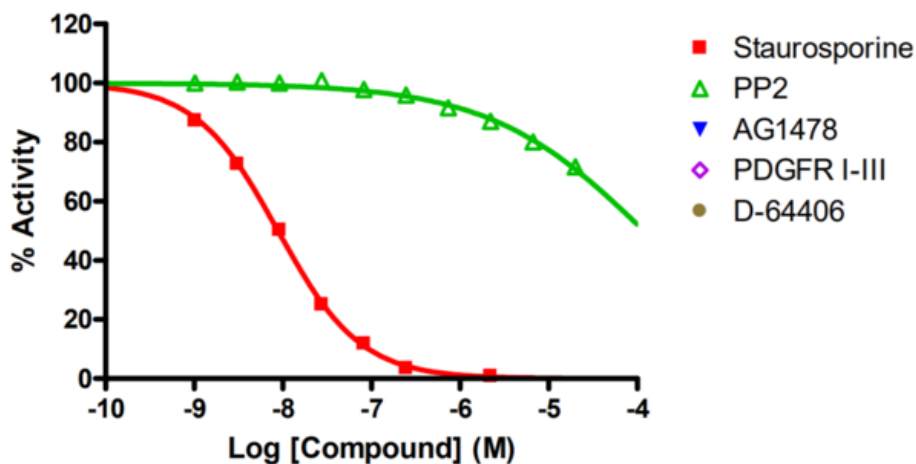
Target Overview

CAT#	IGF1R
Enzyme	Human IGF1R
HGNC	IGF1R
Synonym	JTK13
Substrates	Peptide substrate, [KKKSPGEYVNIEFG], 20 μM ATP 10 μM

Assay Properties

Reaction	Substrate + [³² P]-ATP = ³² P-Substrate + ADP						
Readout	Enzymatic activity						
Reference Compound IC50s	<table border="1"> <thead> <tr> <th>Compounds</th> <th>IC50 (nM)</th> </tr> </thead> <tbody> <tr> <td>Staurosporine</td> <td>9.0</td> </tr> <tr> <td>PP2</td> <td>>20,000</td> </tr> </tbody> </table>	Compounds	IC50 (nM)	Staurosporine	9.0	PP2	>20,000
Compounds	IC50 (nM)						
Staurosporine	9.0						
PP2	>20,000						
Scientific Information	P08069 (http://www.uniprot.org/uniprot/P08069)						
Screening Location	Malvern, USA						

Reference compound IC50 for IGF1R



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/igf1r_kin_malvern

