

DDR2 (T654M) Kinase Assay Service

Activity-based DDR2 (T654M) kinase assay for compound screening and profiling via [radiometric HotSpot™ kinase assay](#).

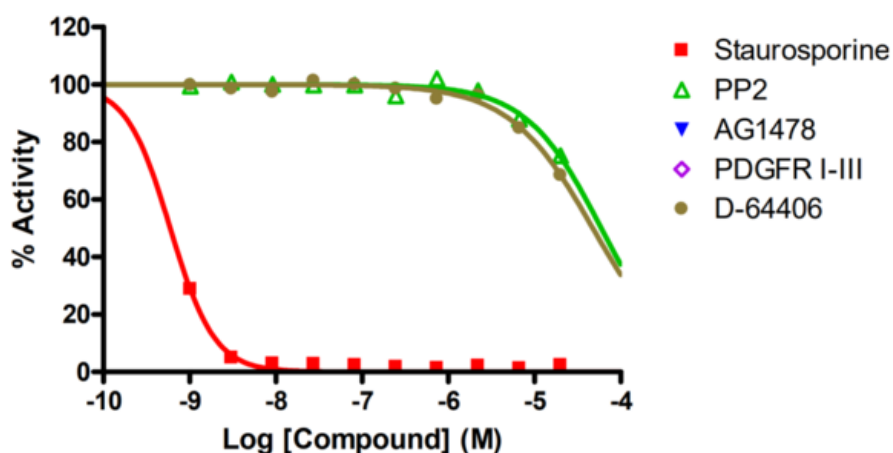
Target Overview

CAT#	DDR2 (T654M)
Enzyme	Human DDR2 (T654M)
HGNC	DDR2 (T654M)
Synonym	MIG20a, NTRKR3, TKT, TYRO10
Substrates	Peptide substrate, [KKSRGDYMTMQIG], 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>0.6</td> </tr> <tr> <td>PP2</td> <td>60,000</td> </tr> <tr> <td>AG1478</td> <td>>100,000</td> </tr> <tr> <td>PDGFR I-III</td> <td>>100,000</td> </tr> <tr> <td>D-64406</td> <td>48,000</td> </tr> </tbody> </table>	Compounds	IC50 (nM)	Staurosporine	0.6	PP2	60,000	AG1478	>100,000	PDGFR I-III	>100,000	D-64406	48,000
Compounds	IC50 (nM)												
Staurosporine	0.6												
PP2	60,000												
AG1478	>100,000												
PDGFR I-III	>100,000												
D-64406	48,000												
Scientific Information	Q16832 (http://www.uniprot.org/uniprot/Q16832)												
Screening Location	Malvern, USA												

Reference compound IC50 for DDR2 (T654M)



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/ddr2_t654m_kin_malvern

