

TIE2 (A1124V) Kinase Assay Service

Activity-based TIE2 (A1124V) kinase assay for compound screening and profiling via [radiometric HotSpot™ kinase assay](#).

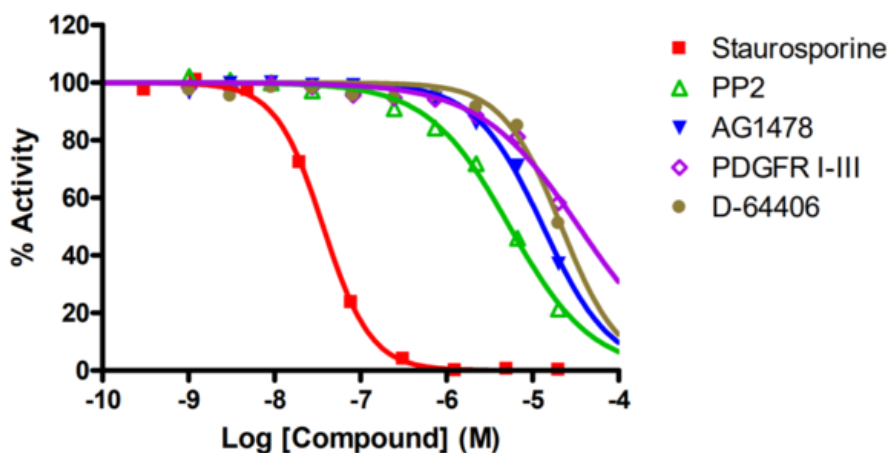
Target Overview

CAT#	TIE2 (A1124V)
Enzyme	Human TIE2 (A1124V)
HGNC	TEK (A1124V)
Synonym	VMCM, TIE-2, VMCM1, CD202B
Substrates	Peptide substrate, poly[Glu:Tyr] (4:1), 0.2 mg/ml 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>37</td> </tr> <tr> <td>PP2</td> <td>5,400</td> </tr> <tr> <td>AG1478</td> <td>13,000</td> </tr> <tr> <td>PDGFR I-III</td> <td>35,000</td> </tr> <tr> <td>D-64406</td> <td>21,000</td> </tr> </tbody> </table>	Compounds	IC50 (nM)	Staurosporine	37	PP2	5,400	AG1478	13,000	PDGFR I-III	35,000	D-64406	21,000
Compounds	IC50 (nM)												
Staurosporine	37												
PP2	5,400												
AG1478	13,000												
PDGFR I-III	35,000												
D-64406	21,000												
Scientific Information	Q02763 (http://www.uniprot.org/uniprot/Q02763)												
Screening Location	Malvern, USA												

Reference compound IC50 for TIE2 (A1124V)



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/tie2_a1124v_kin_malvern

