

c-KIT (V559D) Kinase Assay Service

Activity-based c-KIT (V559D) kinase assay for compound screening and profiling via [radiometric HotSpot™ kinase assay](#).

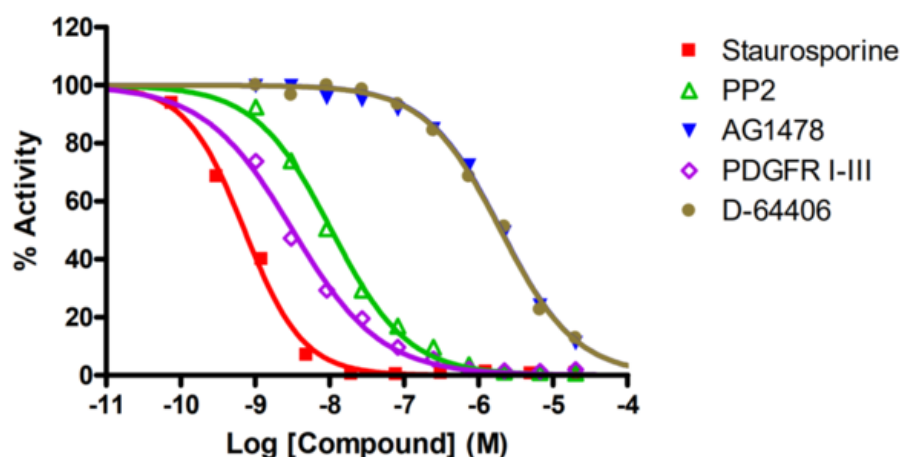
Target Overview

CAT#	c-Kit (V559D)
Enzyme	Human c-Kit (V559D)
HGNC	KIT (V559D)
Synonym	C-Kit, PBT, CD117, SCFR,
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>0.7</td> </tr> <tr> <td>PP2</td> <td>11</td> </tr> <tr> <td>AG1478</td> <td>2,000</td> </tr> <tr> <td>PDGFR I-III</td> <td>3.2</td> </tr> <tr> <td>D-64406</td> <td>2,000</td> </tr> </tbody> </table>	Compounds	IC50 (nM)	Staurosporine	0.7	PP2	11	AG1478	2,000	PDGFR I-III	3.2	D-64406	2,000
Compounds	IC50 (nM)												
Staurosporine	0.7												
PP2	11												
AG1478	2,000												
PDGFR I-III	3.2												
D-64406	2,000												
Scientific Information	P10721 (http://www.uniprot.org/uniprot/P10721)												
Screening Location	Malvern, USA												

Reference compound IC50 for c-Kit (V559D)



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/c-kit_v559d_kin_malvern

