

FGFR3 (K650M) Kinase Assay Service

Activity-based FGFR3 (K650M) kinase assay for compound screening and profiling via [radiometric HotSpot™ kinase assay](#).

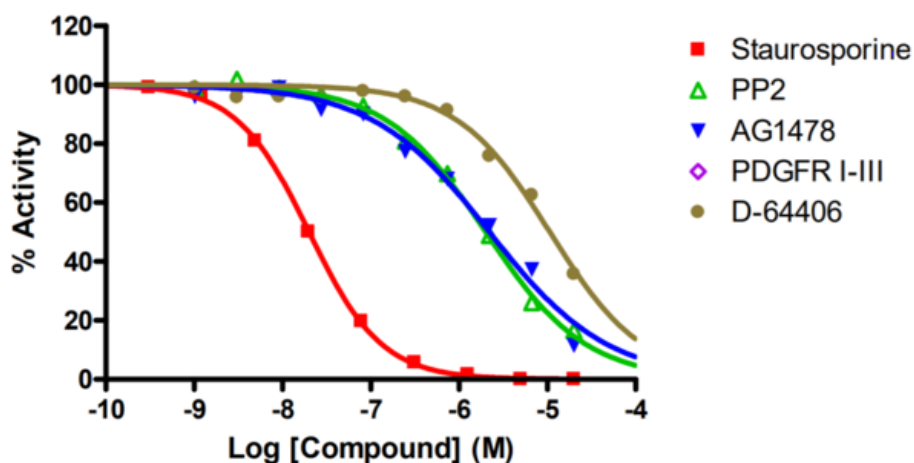
Target Overview

CAT#	FGFR3 (K650M)
Enzyme	Human FGFR3 (K650M)
HGNC	FGFR3 (K650M)
Synonym	ACH, CD333, CEK2, HSFGR3EX, JTK4
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>20</td> </tr> <tr> <td>PP2</td> <td>2,000</td> </tr> <tr> <td>AG1478</td> <td>2,200</td> </tr> <tr> <td>PDGFR I-III</td> <td>>50,000</td> </tr> <tr> <td>D-64406</td> <td>11,000</td> </tr> </tbody> </table>	Compounds	IC50 (nM)	Staurosporine	20	PP2	2,000	AG1478	2,200	PDGFR I-III	>50,000	D-64406	11,000
Compounds	IC50 (nM)												
Staurosporine	20												
PP2	2,000												
AG1478	2,200												
PDGFR I-III	>50,000												
D-64406	11,000												
Scientific Information	P22607 (http://www.uniprot.org/uniprot/P22607)												
Screening Location	Malvern, USA												

Reference compound IC50 for FGFR3 (K650M)



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/fgfr3_k650m_kin_malvern

