

FLT3 (Y591_V592insVDFREYEYD) Kinase Assay Service

Activity-based FLT3 (Y591_V592insVDFREYEYD) kinase assay for compound screening and profiling via [radiometric HotSpot™ kinase assay](#).

Target Overview

CAT#	FLT3 (Y591_V592insVDFREYEYD)
Enzyme	Human FLT3 (Y591_V592insVDFREYEYD)
HGNC	FLT3 (Y591_V592insVDFREYEYD)
Synonym	FLK2, STK1, CD135
Substrates	Peptide substrate, [EAIYAAPFAKKK], 20 μM ATP, 10 μM

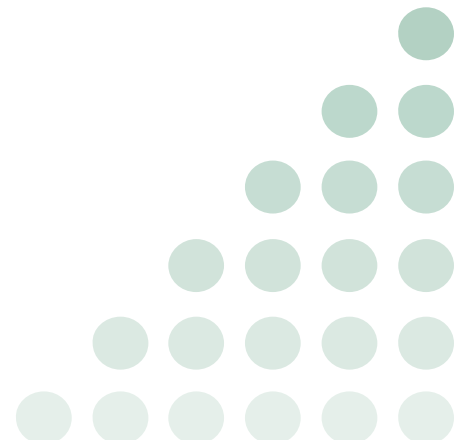
Assay Properties

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>0.9</td> </tr> <tr> <td>PP2</td> <td>2,300</td> </tr> <tr> <td>AG1478</td> <td>190</td> </tr> <tr> <td>PDGFR I-III</td> <td>150</td> </tr> <tr> <td>D-64406</td> <td>18</td> </tr> </tbody> </table>	Compounds	IC50 (nM)	Staurosporine	0.9	PP2	2,300	AG1478	190	PDGFR I-III	150	D-64406	18
Compounds	IC50 (nM)												
Staurosporine	0.9												
PP2	2,300												
AG1478	190												
PDGFR I-III	150												
D-64406	18												
Scientific Information	P36888 (http://www.uniprot.org/uniprot/P36888)												
Screening Location	Malvern, USA												

Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/flt3_y591_v592insvdfreyeyd_kin_malvern



Reference compound IC50 for FLT3 (Y591_V592insVDFREYEYD)

