

ALK (R1275Q) Kinase Assay Service

Activity-based ALK (R1275Q) kinase assay for compound screening and profiling via [radiometric HotSpot™ kinase assay](#).

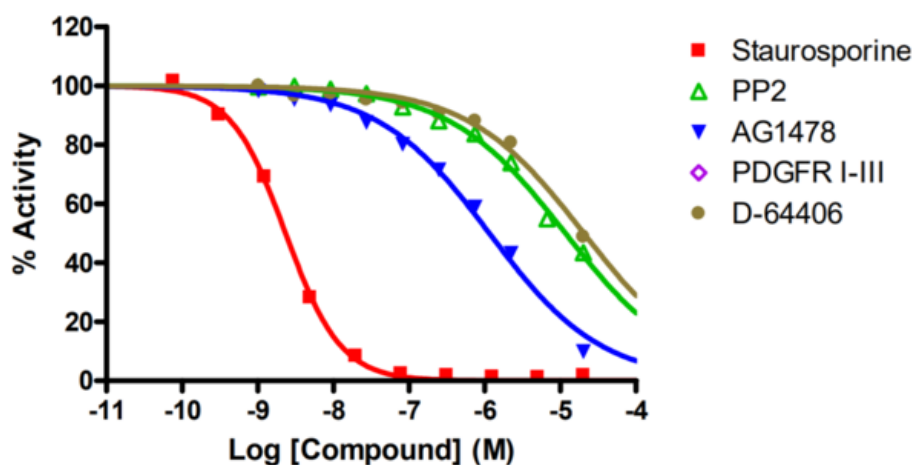
Target Overview

CAT#	ALK (R1275Q)
Enzyme	Human ALK (R1275Q)
HGNC	ALK (R1275Q)
Synonym	CD246, NBLST3
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>2.3</td> </tr> <tr> <td>PP2</td> <td>12,000</td> </tr> <tr> <td>AG1478</td> <td>1,100</td> </tr> <tr> <td>PDGFR I-III</td> <td>>100,000</td> </tr> <tr> <td>D-64406</td> <td>20,000</td> </tr> </tbody> </table>	Compounds	IC50 (nM)	Staurosporine	2.3	PP2	12,000	AG1478	1,100	PDGFR I-III	>100,000	D-64406	20,000
Compounds	IC50 (nM)												
Staurosporine	2.3												
PP2	12,000												
AG1478	1,100												
PDGFR I-III	>100,000												
D-64406	20,000												
Scientific Information	Q9UM73 (http://www.uniprot.org/uniprot/Q9UM73)												
Screening Location	Malvern, USA												

Reference compound IC50 for ALK (R1275Q)



Let's Discover Together.

© Aug 2022 Reaction Biology. All rights reserved.

https://www.reactionbiology.com/datasheet/alk_r1275q_kin_malvern

