

ALK-TPM1 Kinase Assay Service

Activity-based ALK-TPM1 kinase assay for compound screening and profiling via [radiometric HotSpot™ kinase assay](#).

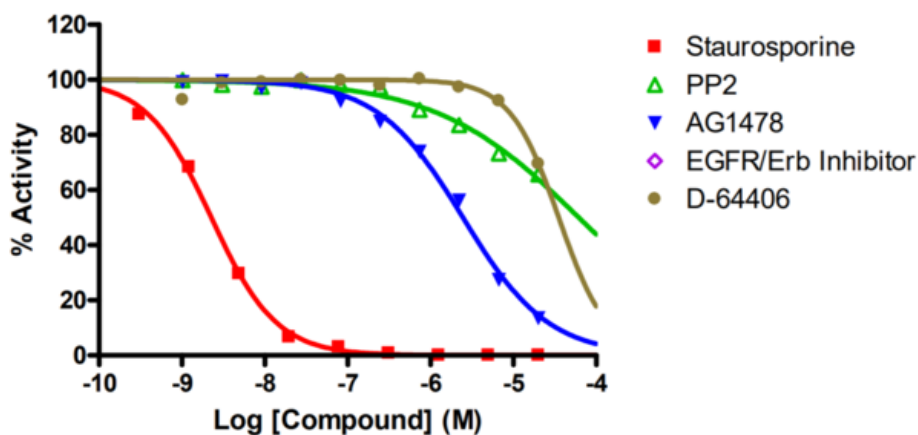
Target Overview

CAT#	ALK, TPM1-
Enzyme	Human TPM1-ALK fusion
HGNC	ALK-TPM1
Synonym	CD246, NBLST3
Substrates	Peptide substrate, poly[Glu:Tyr] (4:1), 0.2 mg/ml ATP 10 μM

Assay Properties

Reaction	Substrate + [³³ P]-ATP = ³³ P-Substrate + ADP	
Readout	Enzymatic activity	
Reference Compound IC50s	Compounds	IC50 (nM)
	Staurosporine	2.3
	PP2	62,000
	AG1478	2,500
	EGFR/Erb Inhibitor	>100,000
	D-64406	35,000
Scientific Information	TPM1; P09493 (http://www.uniprot.org/uniprot/P09493)ALK; Q9UM73 (http://www.uniprot.org/uniprot/Q9UM73)	
Screening Location	Malvern, USA	

Reference compound IC50 for ALK-TPM1



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/alk-tpm1_kin_malvern

