

TRKA (F589L) Kinase Assay Service

Activity-based TRKA (F589L) kinase assay for compound screening and profiling via [radiometric HotSpot™ kinase assay](#).

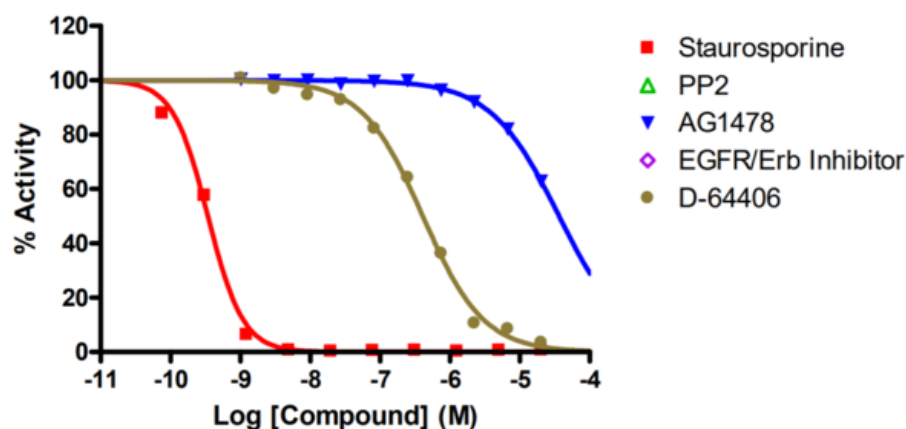
Target Overview

CAT#	TRKA (F589L)
Enzyme	Human TRKA (F589L)
HGNC	NTRK1 (F589L)
Synonym	MTC, TRK, TRK1
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	Compounds	IC50 (nM)
	Staurosporine	0.34
	PP2	>100,000
	AG1478	36,000
	EGFR/Erb Inhibitor	>100,000
	D-64406	410
Scientific Information	P04629 (http://www.uniprot.org/uniprot/P04629)	
Screening Location	Malvern, USA	

Reference compound IC50 for TRKA (F589L)



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/trka_f589l_kin_malvern

