

RET (S904F) Kinase Assay Service

Activity-based RET (S904F) kinase assay for compound screening and profiling via [radiometric HotSpot™ kinase assay](#).

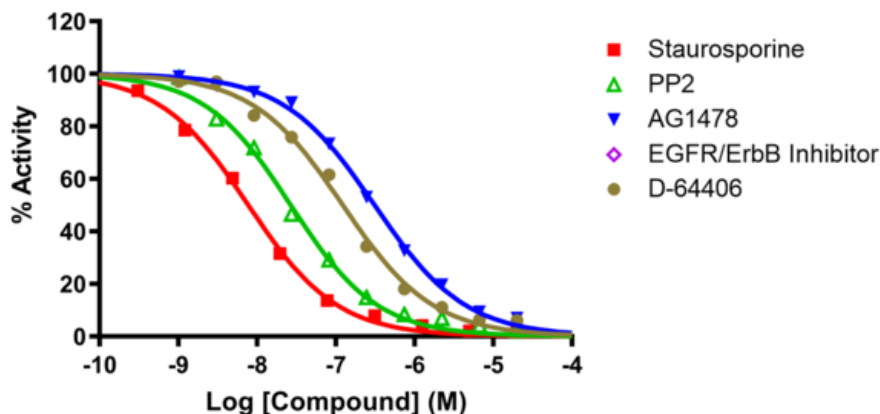
Target Overview

CAT#	RET (S904F)
Enzyme	Human RET (S904F)
HGNC	RET (S904F)
Synonym	CDHF12, CDHR16, HSCR1, MEN2A, MEN2B, MTC1, PTC, RET51
Substrates	Peptide substrate, [RRRRRRRRRRRVYSTDYRLFNPS], 20 μM ATP 10 μM

Assay Properties

Reaction	Substrate + [³² P]-ATP = ³² P-Substrate + ADP	
Readout	Enzymatic activity	
Reference Compound IC50s	Compounds	IC50 (nM)
	Staurosporine	7.7
	PP2	27
	AG1478	310
	EGFR/ErbB Inhibitor	>100,000
	D-64406	120
Scientific Information	P07949 (http://www.uniprot.org/uniprot/P07949)	
Screening Location	Malvern, USA	

Reference compound IC50 for RET (S904F)



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/ret_s904f_kin_malvern

