

# RET (E762Q) Kinase Assay Service

Activity-based RET (E762Q) kinase assay for compound screening and profiling via [radiometric HotSpot™ kinase assay](#).

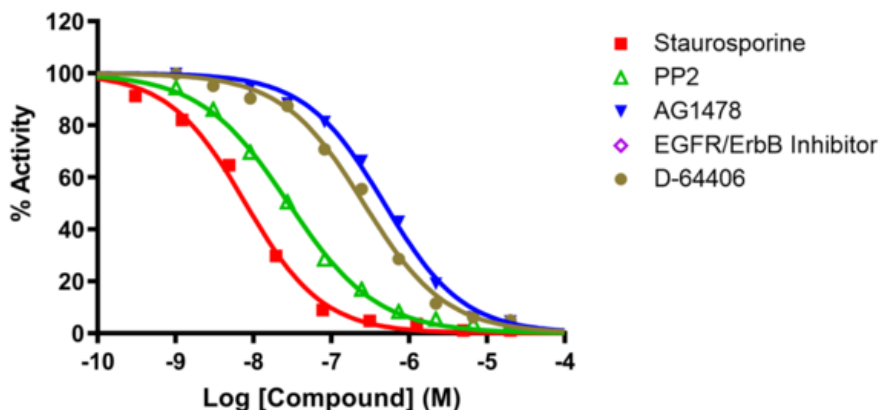
## Target Overview

<b>CAT#</b>	RET (E762Q)
<b>Enzyme</b>	Human RET (E762Q)
<b>HGNC</b>	RET (E762Q)
<b>Synonym</b>	CDHF12, CDHR16, HSCR1, MEN2A, MEN2B, MTC1, PTC, RET51
<b>Substrates</b>	Peptide substrate, [RRRRRRRRRRRVYSTDYRLFNPS], 20 μM ATP 10 μM

## Assay Properties

<b>Reaction</b>	Substrate + [ <sup>32</sup> P]-ATP = <sup>32</sup> P-Substrate + ADP	
<b>Readout</b>	Enzymatic activity	
<b>Reference Compound IC50s</b>	<b>Compounds</b>	<b>IC50 (nM)</b>
	Staurosporine	8.0
	PP2	28
	AG1478	480
	EGFR/ErbB Inhibitor	>100,000
	D-64406	260
<b>Scientific Information</b>	<a href="http://www.uniprot.org/uniprot/P07949">P07949</a> ( <a href="http://www.uniprot.org/uniprot/P07949">http://www.uniprot.org/uniprot/P07949</a> )	
<b>Screening Location</b>	Malvern, USA	

Reference compound IC50 for RET (E762Q)



Let's Discover Together.

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

[https://www.reactionbiology.com/datasheet/ret\\_e762q\\_kin\\_malvern](https://www.reactionbiology.com/datasheet/ret_e762q_kin_malvern)

