

c-MET (H1094Y) Kinase Assay Service

Activity-based c-MET (H1094Y) kinase assay for compound screening and profiling via [radiometric HotSpot™ kinase assay](#).

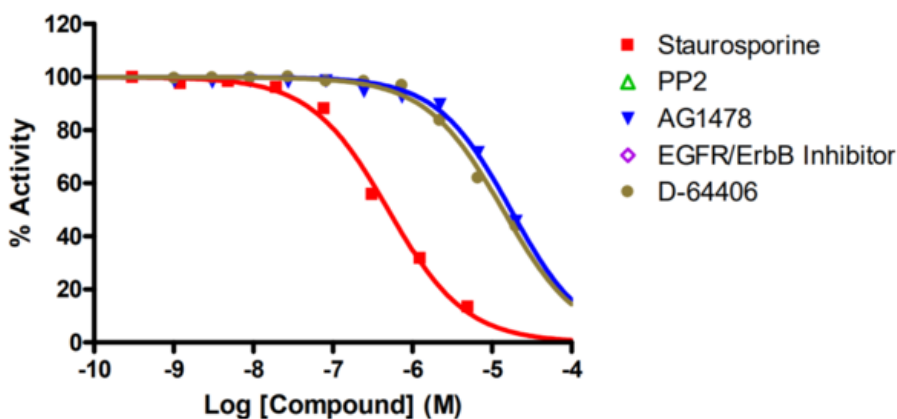
Target Overview

CAT#	c-MET (H1094Y)
Enzyme	Human c-MET (H1094Y)
HGNC	MET (H1094Y)
Synonym	c-MET, HGFR, RCCP2
Substrates	MBP (Myelin Basic Protein), 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	500
	PP2	>100,000
	AG1478	17,000
	EGFR/Erb Inhibitor	>100,000
	D-64406	14,000
Scientific Information	P08581 (http://www.uniprot.org/uniprot/P08581)	
Screening Location	Malvern, USA	

Reference compound IC50 for C-MET (H1094Y)



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/c-met_h1094y_kin_malvern

