

EGFR (d746-750/T790M/C797S/L858R) Kinase Assay Service

Activity-based EGFR (d746-750/T790M/C797S/L858R) kinase assay for compound screening and profiling via [radiometric HotSpot™ kinase assay](#).

Target Overview

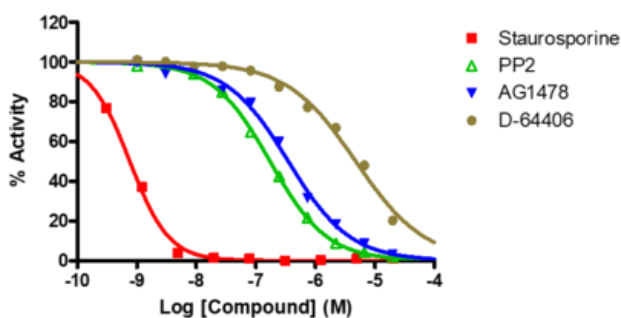
CAT#	EGFR (d746-750/T790M/C797S/L858R)
Enzyme	Human EGFR (d746-750/ T790M/C797S/L858R)
HGNC	EGFR (d746-750/T790M/C797S/L858R)
Synonym	Epidermal growth factor receptor, Receptor tyrosine-protein kinase ErbB-1, HER1, EGF-R, ERBB, mENA, ERBB1
Substrates	Peptide substrate, poly[Glu:Tyr] (4:1), 0.2 mg/ml ATP 10 μM

Assay Properties

Reaction	Substrate + [³³ P]-ATP = ³³ P-Substrate + ADP	
Readout	Enzymatic activity	
Reference Compound IC50s	Compounds	IC50 (nM)
	Staurosporine	0.8
	PP2	174
	AG1478	350
	EGFR/ErbB Inhibitor	>100,000
	D-64406	4,800
Scientific Information	P00533 (http://www.uniprot.org/uniprot/P00533)	
Screening Location	Malvern, USA	

Reference compound IC50 for EGFR (d746-750/T790M/C797S/L858R)

Reference compound IC50 for EGFR (d746-750/T790M/C797S/L858R)



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/egfr_d746-750_t790m_c797s_l858r_kin_malvern

