

EGFR (d746-750/C797S) Kinase Assay Service

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

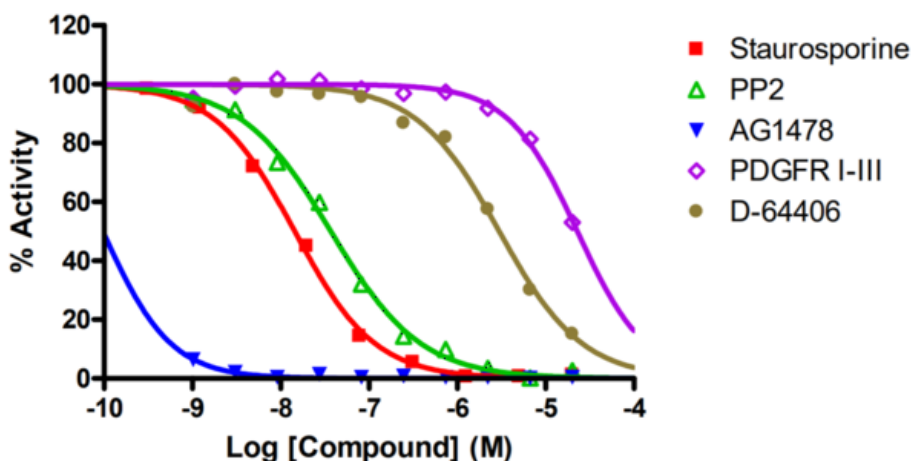
Target Overview

CAT#	EGFR (d746-750/C797S)
Enzyme	Human EGFR (d746-750/C797S)
HGNC	EGFR (d746-750/C797S)
Synonym	ERBB, mENA, ERBB1
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	<table border="1"> <thead> <tr> <th>Compounds</th> <th>IC50 (nM)</th> </tr> </thead> <tbody> <tr> <td>Staurosporine</td> <td>14</td> </tr> <tr> <td>PP2</td> <td>37</td> </tr> <tr> <td>AG1478</td> <td><1.0</td> </tr> <tr> <td>PDGFR I-III</td> <td>22,800</td> </tr> <tr> <td>D-64406</td> <td>3,000</td> </tr> </tbody> </table>	Compounds	IC50 (nM)	Staurosporine	14	PP2	37	AG1478	<1.0	PDGFR I-III	22,800	D-64406	3,000
Compounds	IC50 (nM)												
Staurosporine	14												
PP2	37												
AG1478	<1.0												
PDGFR I-III	22,800												
D-64406	3,000												
Scientific Information	P00533 (http://www.uniprot.org/uniprot/P00533)												
Screening Location	Malvern, USA												

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



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/egfr_d746-750_c797s_kin_malvern

