

EGFR Kinase Assay Service

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

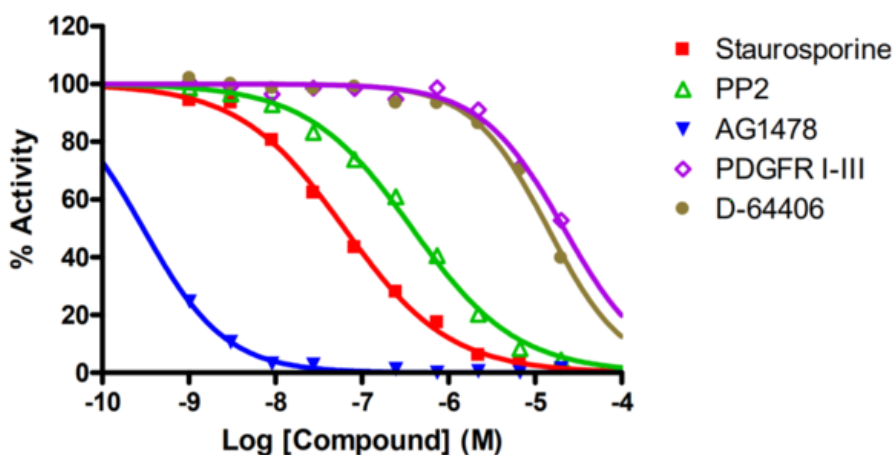
Target Overview

CAT#	EGFR
Enzyme	Human EGFR
HGNC	EGFR
Synonym	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	<table border="1"> <thead> <tr> <th>Compounds</th> <th>IC50 (nM)</th> </tr> </thead> <tbody> <tr> <td>Staurosporine</td> <td>63</td> </tr> <tr> <td>PP2</td> <td>380</td> </tr> <tr> <td>AG1478</td> <td><0.1</td> </tr> <tr> <td>PDGFR I-III</td> <td>21,000</td> </tr> <tr> <td>D-64406</td> <td>14,000</td> </tr> </tbody> </table>	Compounds	IC50 (nM)	Staurosporine	63	PP2	380	AG1478	<0.1	PDGFR I-III	21,000	D-64406	14,000
Compounds	IC50 (nM)												
Staurosporine	63												
PP2	380												
AG1478	<0.1												
PDGFR I-III	21,000												
D-64406	14,000												
Scientific Information	P00533 (http://www.uniprot.org/uniprot/P00533)												
Screening Location	Malvern, USA												

Reference compound IC50 for EGFR



Let's Discover Together.

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

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

