

# EGFR (K716A/C797S/L858R) Kinase Assay Service

Activity-based EGFR (K716A/C797S/L858R) kinase assay for compound screening and profiling via [radiometric HotSpot™ kinase assay](#).

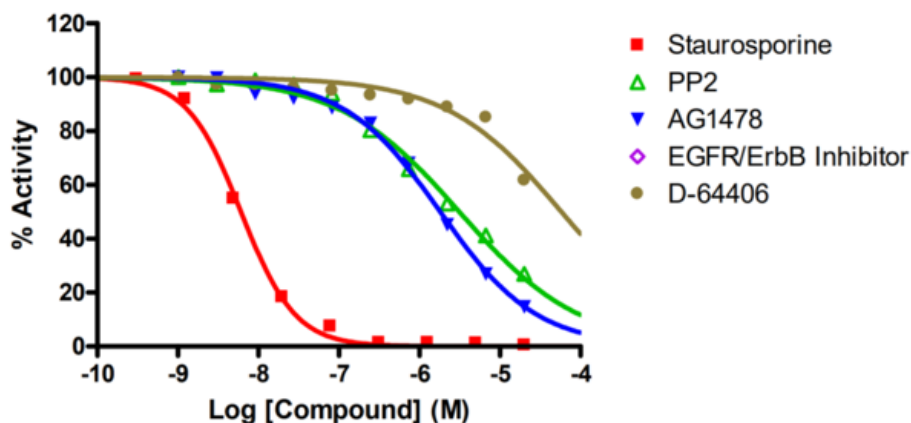
## Target Overview

<b>CAT#</b>	EGFR (K716A/C797S/L858R)
<b>Enzyme</b>	Human EGFR (K716A/C797S/L858R)
<b>HGNC</b>	EGFR (K716A/C797S/L858R)
<b>Synonym</b>	ERBB, mENA, ERBB1
<b>Substrates</b>	Peptide substrate, poly[Glu:Tyr] (4:1), 0.2 mg/ml 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	6.1
	PP2	3,100
	AG1478	1,800
	EGFR/ErbB Inhibitor	>100,000
	D-64406	58,000
<b>Scientific Information</b>	<a href="http://www.uniprot.org/uniprot/P00533">P00533</a> ( <a href="http://www.uniprot.org/uniprot/P00533">http://www.uniprot.org/uniprot/P00533</a> )	
<b>Screening Location</b>	Malvern, USA	

## Reference compound IC50 for EGFR (K716A/C797S/L858R)



Let's Discover Together.

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

[https://www.reactionbiology.com/datasheet/egfr\\_k716a\\_c797s\\_l858r\\_kin\\_malvern](https://www.reactionbiology.com/datasheet/egfr_k716a_c797s_l858r_kin_malvern)

