

EPHB3 Kinase Assay Service

Activity-based EPHB3 kinase assay for compound screening and profiling via [radiometric HotSpot™ kinase assay](#).

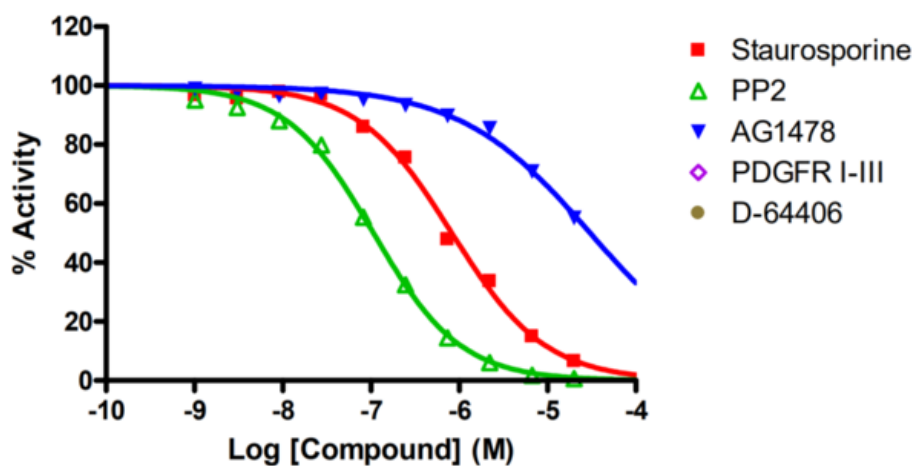
Target Overview

CAT#	EPHB3
Enzyme	Human EPHB3
HGNC	EPHB3
Synonym	ETK2, HEK2, TYRO6
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>825</td> </tr> <tr> <td>PP2</td> <td>107</td> </tr> <tr> <td>AG1478</td> <td>30,000</td> </tr> </tbody> </table>	Compounds	IC50 (nM)	Staurosporine	825	PP2	107	AG1478	30,000
Compounds	IC50 (nM)								
Staurosporine	825								
PP2	107								
AG1478	30,000								
Scientific Information	P54753 (http://www.uniprot.org/uniprot/P54753)								
Screening Location	Malvern, USA								

Reference compound IC50 for EPHB3



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/ephb3_kin_malvern

