

BMX/ETK Kinase Assay Service

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

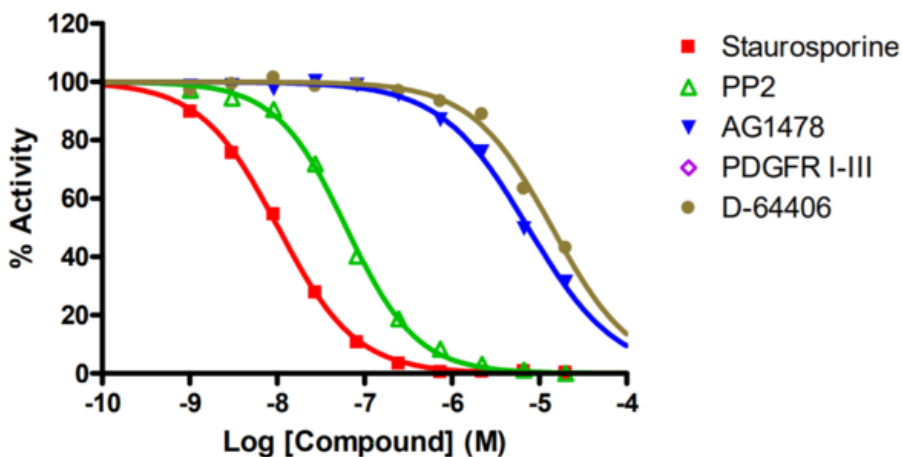
Target Overview

CAT#	BMX
Enzyme	Human BMX
HGNC	BMX
Synonym	ETK, PSCTK2, PSCTK3
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>10</td> </tr> <tr> <td>PP2</td> <td>62</td> </tr> <tr> <td>AG1478</td> <td>7,400</td> </tr> <tr> <td>D-64406</td> <td>14,000</td> </tr> </tbody> </table>	Compounds	IC50 (nM)	Staurosporine	10	PP2	62	AG1478	7,400	D-64406	14,000
Compounds	IC50 (nM)										
Staurosporine	10										
PP2	62										
AG1478	7,400										
D-64406	14,000										
Scientific Information	P51813 (http://www.uniprot.org/uniprot/P51813)										
Screening Location	Malvern, USA										

Reference compound IC50 for BMX/ETK



Let's Discover Together.

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

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

