

BLK Kinase Assay Service

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

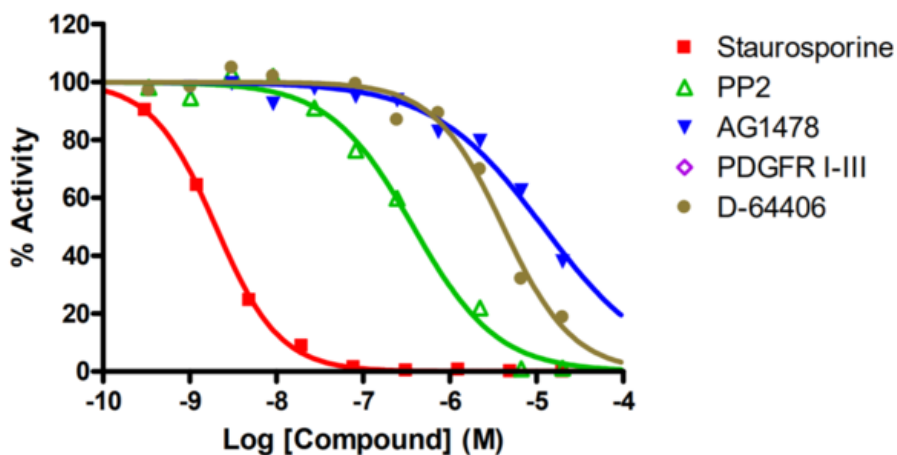
Target Overview

CAT#	BLK
Enzyme	Human BLK
HGNC	BLK
Synonym	BLK
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>2.1</td> </tr> <tr> <td>PP2</td> <td>379</td> </tr> <tr> <td>AG1478</td> <td>12,000</td> </tr> <tr> <td>D-64406</td> <td>4,100</td> </tr> </tbody> </table>	Compounds	IC50 (nM)	Staurosporine	2.1	PP2	379	AG1478	12,000	D-64406	4,100
Compounds	IC50 (nM)										
Staurosporine	2.1										
PP2	379										
AG1478	12,000										
D-64406	4,100										
Scientific Information	P51451 (http://www.uniprot.org/uniprot/P51451)										
Screening Location	Malvern, USA										

Reference compound IC50 for BLK



Let's Discover Together.

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

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

