

PRKX Kinase Assay Service

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

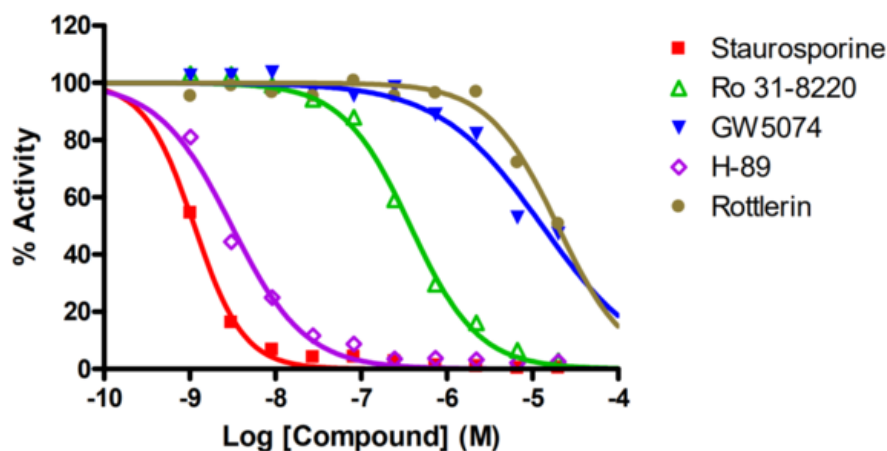
Target Overview

CAT#	PRKX
Enzyme	Human PRKX
HGNC	PRKX
Synonym	PKX1
Substrates	Peptide substrate, [LRRASLG], 100 μM 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>1.1</td> </tr> <tr> <td>Ro 31-8220</td> <td>380</td> </tr> <tr> <td>GW5074</td> <td>13,000</td> </tr> <tr> <td>H-89</td> <td>3.1</td> </tr> <tr> <td>Rottlerin</td> <td>19,700</td> </tr> </tbody> </table>	Compounds	IC50 (nM)	Staurosporine	1.1	Ro 31-8220	380	GW5074	13,000	H-89	3.1	Rottlerin	19,700
Compounds	IC50 (nM)												
Staurosporine	1.1												
Ro 31-8220	380												
GW5074	13,000												
H-89	3.1												
Rottlerin	19,700												
Scientific Information	P51817 (http://www.uniprot.org/uniprot/P51817)												
Screening Location	Malvern, USA												

Reference compound IC50 for PRKX



Let's Discover Together.

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

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

