

PKD2/PRKD2 Kinase Assay Service

Activity-based PKD2/PRKD2 kinase assay for compound screening and profiling via [radiometric HotSpot™ kinase assay](#).

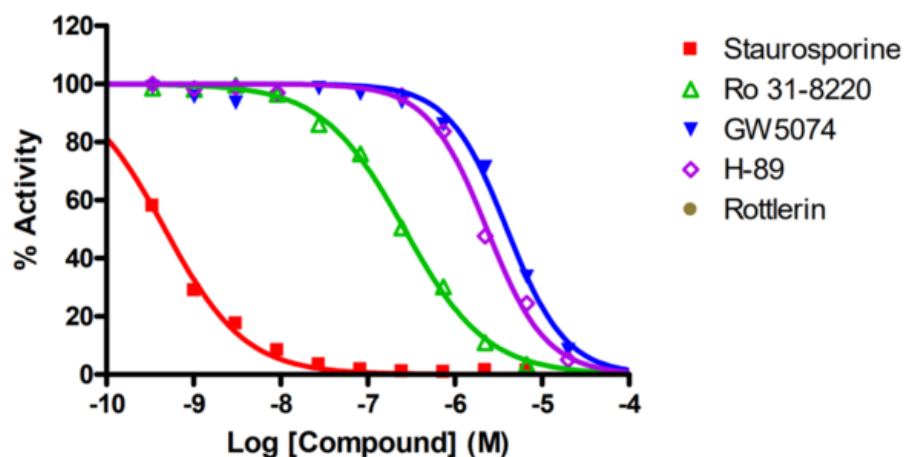
Target Overview

CAT#	PKD2
Enzyme	Human PKD2
HGNC	PRKD2
Synonym	PKD2, HSPC187
Substrates	Peptide substrate, [KKLNRTLSVA], 20 μM 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>0.47</td> </tr> <tr> <td>Ro 31-8220</td> <td>263</td> </tr> <tr> <td>GW 5074</td> <td>4,000</td> </tr> <tr> <td>H-89</td> <td>2,300</td> </tr> </tbody> </table>	Compounds	IC50 (nM)	Staurosporine	0.47	Ro 31-8220	263	GW 5074	4,000	H-89	2,300
Compounds	IC50 (nM)										
Staurosporine	0.47										
Ro 31-8220	263										
GW 5074	4,000										
H-89	2,300										
Scientific Information	Q9BZL6 (http://www.uniprot.org/uniprot/Q9BZL6)										
Screening Location	Malvern, USA										

Reference compound IC50 for PKD2/PRKD2



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/pkd2_prkd2_kin_malvern

