

PKG1a Kinase Assay Service

Activity-based PKG1a kinase assay for compound screening and profiling via [radiometric HotSpot™ kinase assay](#).

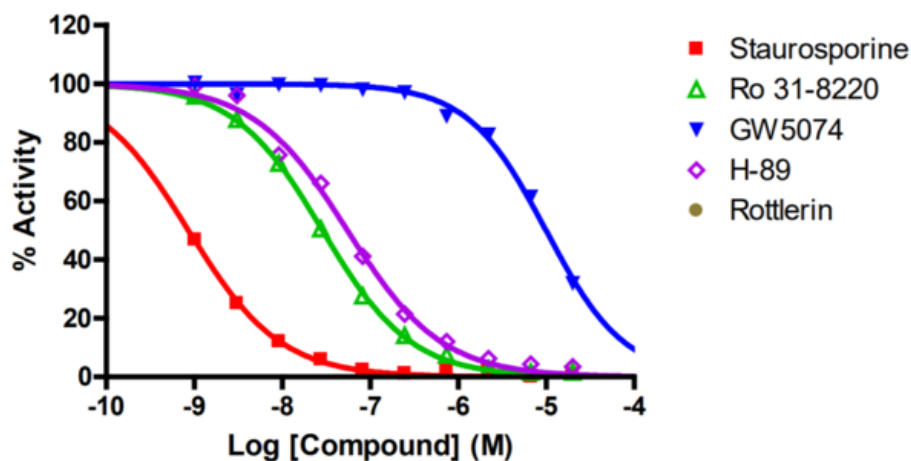
Target Overview

CAT#	PKG1
Enzyme	Human PKG1
HGNC	PRKG1
Synonym	PKG1, PGK, PRKG1B, PRKGR1B
Substrates	Peptide substrate, [LRRASLG], 100 μM ATP 1 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>0.86</td> </tr> <tr> <td>Ro 31-8220</td> <td>29</td> </tr> <tr> <td>GW 5074</td> <td>10,000</td> </tr> <tr> <td>H-89</td> <td>55</td> </tr> </tbody> </table>	Compounds	IC50 (nM)	Staurosporine	0.86	Ro 31-8220	29	GW 5074	10,000	H-89	55
Compounds	IC50 (nM)										
Staurosporine	0.86										
Ro 31-8220	29										
GW 5074	10,000										
H-89	55										
Scientific Information	Q13976 (http://www.uniprot.org/uniprot/Q13976)										
Screening Location	Malvern, USA										

Reference compound IC50 for PKG1



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/pkg1a_kin_malvern

