

VRK1 Kinase Assay Service

Activity-based VRK1 kinase assay for compound screening and profiling via [radiometric HotSpot™ kinase assay](#).

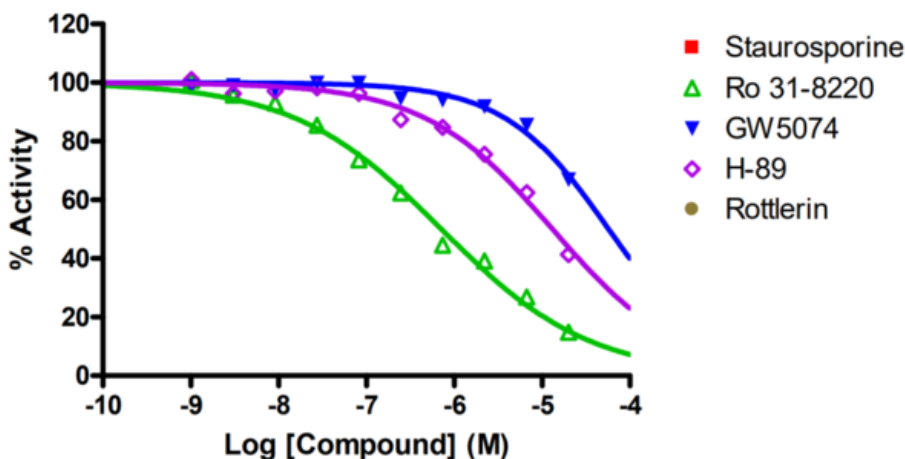
Target Overview

CAT#	VRK1
Enzyme	Human VRK1
HGNC	VRK1
Synonym	VRK1
Substrates	Peptide substrate, [KEAKEKRQEIQAKRRRLSSLRASTSKSGGSQK], 20 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>>100,000</td> </tr> <tr> <td>Ro 31-8220</td> <td>720</td> </tr> <tr> <td>GW 5074</td> <td>57,000</td> </tr> <tr> <td>H-89</td> <td>13,000</td> </tr> <tr> <td>Rottlerin</td> <td>>100,000</td> </tr> </tbody> </table>	Compounds	IC50 (nM)	Staurosporine	>100,000	Ro 31-8220	720	GW 5074	57,000	H-89	13,000	Rottlerin	>100,000
Compounds	IC50 (nM)												
Staurosporine	>100,000												
Ro 31-8220	720												
GW 5074	57,000												
H-89	13,000												
Rottlerin	>100,000												
Scientific Information	Q99986 (http://www.uniprot.org/uniprot/Q99986)												
Screening Location	Malvern, USA												

Reference compound IC50 for VRK1



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/vrk1_kin_malvern

