

ROCK1 Kinase Assay Service

Activity-based ROCK1 kinase assay for compound screening and profiling via [radiometric HotSpot™ kinase assay](#).

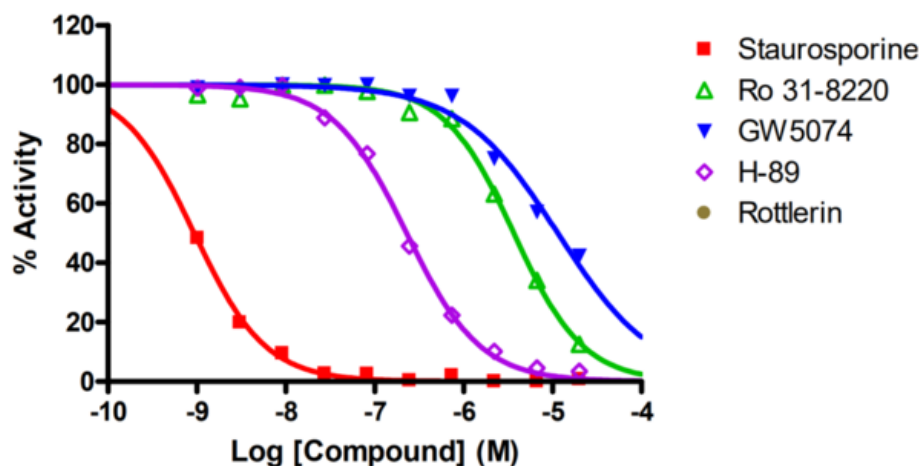
Target Overview

CAT#	ROCK1
Enzyme	Human ROCK1
HGNC	ROCK1
Synonym	P160ROCK
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>0.93</td> </tr> <tr> <td>Ro 31-8220</td> <td>3,700</td> </tr> <tr> <td>GW 5074</td> <td>11,500</td> </tr> <tr> <td>H-89</td> <td>230</td> </tr> </tbody> </table>	Compounds	IC50 (nM)	Staurosporine	0.93	Ro 31-8220	3,700	GW 5074	11,500	H-89	230
Compounds	IC50 (nM)										
Staurosporine	0.93										
Ro 31-8220	3,700										
GW 5074	11,500										
H-89	230										
Scientific Information	Q13464 (http://www.uniprot.org/uniprot/Q13464)										
Screening Location	Malvern, USA										

Reference compound IC50 for ROCK1



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/rock1_kin_malvern

