

SIK1 Kinase Assay Service

Activity-based SIK1 kinase assay for compound screening and profiling via [radiometric HotSpot™ kinase assay](#).

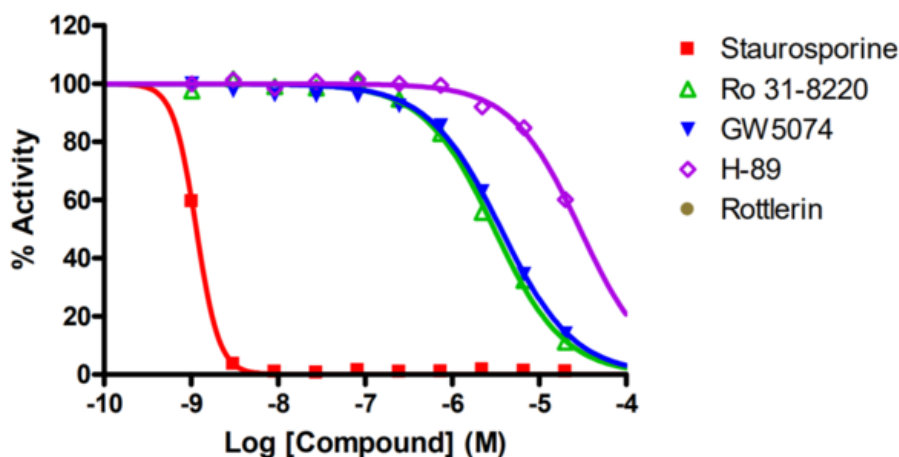
Target Overview

CAT#	SIK1
Enzyme	Human SIK1
HGNC	SIK1
Synonym	SIK, MSK, SNF1LK
Substrates	Peptide substrate, [AMARAASAAALARRR], 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>1.2</td> </tr> <tr> <td>Ro 31-8220</td> <td>3,000</td> </tr> <tr> <td>GW5074</td> <td>3,600</td> </tr> <tr> <td>H-89</td> <td>29,000</td> </tr> <tr> <td>Rottlerin</td> <td>>30,000</td> </tr> </tbody> </table>	Compounds	IC50 (nM)	Staurosporine	1.2	Ro 31-8220	3,000	GW5074	3,600	H-89	29,000	Rottlerin	>30,000
Compounds	IC50 (nM)												
Staurosporine	1.2												
Ro 31-8220	3,000												
GW5074	3,600												
H-89	29,000												
Rottlerin	>30,000												
Scientific Information	P57059 (http://www.uniprot.org/uniprot/P57059)												
Screening Location	Malvern, USA												

Reference compound IC50 for SIK1



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/sik1_kin_malvern

