

# SGK1 Kinase Assay Service

Activity-based SGK1 kinase assay for compound screening and profiling via [radiometric HotSpot™ kinase assay](#).

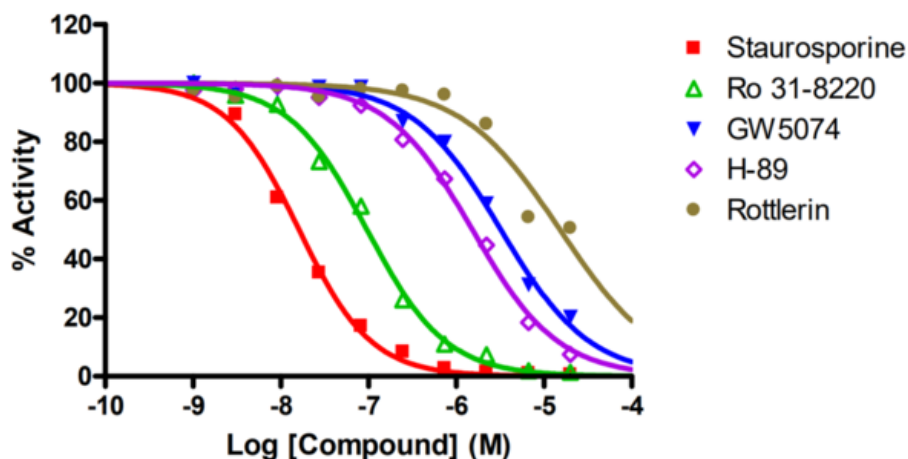
## Target Overview

CAT#	SGK1
Enzyme	Human SGK1
HGNC	SGK1
Synonym	SGK
Substrates	Peptide substrate, [KGSVSGRPRVSSFAEG], 20 μM ATP 10 μM

## Assay Properties

Reaction	Substrate + [ <sup>32</sup> P]-ATP = <sup>32</sup> 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>16</td> </tr> <tr> <td>Ro 31-8220</td> <td>96</td> </tr> <tr> <td>GW 5074</td> <td>3,200</td> </tr> <tr> <td>H-89</td> <td>1,500</td> </tr> <tr> <td>Rottlerin</td> <td>15,000</td> </tr> </tbody> </table>	Compounds	IC50 (nM)	Staurosporine	16	Ro 31-8220	96	GW 5074	3,200	H-89	1,500	Rottlerin	15,000
Compounds	IC50 (nM)												
Staurosporine	16												
Ro 31-8220	96												
GW 5074	3,200												
H-89	1,500												
Rottlerin	15,000												
Scientific Information	<a href="http://www.uniprot.org/uniprot/O00141">O00141</a> ( <a href="http://www.uniprot.org/uniprot/O00141">http://www.uniprot.org/uniprot/O00141</a> )												
Screening Location	Malvern, USA												

## Reference compound IC50 for SGK1



Let's Discover Together.

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

[https://www.reactionbiology.com/datasheet/sgk1\\_kin\\_malvern](https://www.reactionbiology.com/datasheet/sgk1_kin_malvern)

